

PRODUCT INFORMATION



Synthetic Cannabinoid Screening Library (96-Well)

Item No. 9002891 • Batch No. 0698808

Panels are routinely re-evaluated to include new catalog introductions as the research evolves.

Page 1 of 19

Plate	Well	Contents	Item Number
1	A1	Unused	
1	A2	JWH 007	10266
1	A3	MDA 19	10563
1	A4	JWH 251	10578
1	A5	JWH 073 2-hydroxyindole metabolite	10633
1	A6	MDA 77	10639
1	A7	JWH 182	10643
1	A8	JWH 210	10644
1	A9	JWH 098	10680
1	A10	JWH 018 N-(3-methylbutyl) isomer	10691
1	A11	JWH 018 6-methoxyindole analog	10697
1	A12	Unused	
1	B1	Unused	
1	B2	JWH 201	10721
1	B3	JWH 302	10722
1	B4	JWH 200 6-hydroxyindole metabolite	10746
1	B5	(±)-JWH 073 N-(3-hydroxybutyl) metabolite	10795
1	B6	(±)-JWH 018 N-(4-hydroxypentyl) metabolite	10796
1	B7	JWH 307	10797
1	B8	RCS-4-C4 homolog	10798
1	B9	JWH 031	10824
1	B10	JWH 145	10825
1	B11	JWH 370	10827
1	B12	Unused	
1	C1	Unused	
1	C2	JWH 369	10828
1	C3	JWH 368	10829
1	C4	JWH 309	10830
1	C5	JWH 030	10831
1	C6	JWH 073 2-methylnaphthyl analog	10848
1	C7	JWH 016	10849
1	C8	JWH 020	10850
1	C9	AM2201 2'-naphthyl isomer	10862
1	C10	RCS-8 4-methoxy isomer	10863
1	C11	RCS-8 3-methoxy isomer	10864
1	C12	Unused	

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/09/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 • USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 2 of 19

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	RCS-4 2-methoxy isomer	10865
1	D3	RCS-4 3-methoxy isomer	10866
1	D4	JWH 203 4-chlorophenyl isomer	10867
1	D5	JWH 203 3-chlorophenyl isomer	10868
1	D6	AM694 4-iodo isomer	10869
1	D7	AM694 3-iodo isomer	10870
1	D8	JWH 122 N-(5-hydroxypentyl) metabolite	10925
1	D9	RCS-4 N-(5-hydroxypentyl) metabolite	10935
1	D10	RCS-4 N-(4-hydroxypentyl) metabolite	10936
1	D11	RCS-4 N-(5-carboxypentyl) metabolite	10937
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	JWH 250 N-pentanoic acid metabolite	10938
1	E3	JWH 250 N-(4-hydroxypentyl) metabolite	10939
1	E4	JWH 210 N-(4-hydroxypentyl) metabolite	10940
1	E5	JWH 210 N-pentanoic acid metabolite	10941
1	E6	JWH 398 N-(4-hydroxypentyl) metabolite	10943
1	E7	AM2233	11008
1	E8	JWH 018 N-(5-bromopentyl) analog	11047
1	E9	(±)-JWH 018 N-(2-hydroxypentyl) metabolite	11119
1	E10	JWH 249	11153
1	E11	A-834735	11163
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	AM2201 6-hydroxyindole metabolite	11192
1	F3	AM2201 7-hydroxyindole metabolite	11193
1	F4	AM2201 2-hydroxyindole metabolite	11194
1	F5	JWH 175	11201
1	F6	AM1248	11282
1	F7	AM2232	11503
1	F8	AM679	11504
1	F9	AM1220 azepane isomer	11583
1	F10	JWH 018 N-(1-ethylpropyl) isomer	11585
1	F11	JWH 018 2'-naphthyl-N-(1-ethylpropyl) isomer	11586
1	F12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 3 of 19

Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	A-796260	11606
1	G3	JWH 176	11617
1	G4	Cannabipiperidiethanone	11655
1	G5	JWH 213	11659
1	G6	1'-Naphthoyl Indole	11687
1	G7	XLR11 N-(4-pentenyl) analog	11688
1	G8	AB-005	11766
1	G9	UR-144 N-pentanoic acid metabolite	11773
1	G10	UR-144 N-(4-hydroxypentyl) metabolite	11774
1	G11	UR-144 N-(5-hydroxypentyl) metabolite	11775
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	MAM2201 N-(4-hydroxypentyl) metabolite	11778
1	H3	MAM2201 N-pentanoic acid metabolite	11779
1	H4	JWH 122 N-(4-hydroxypentyl) metabolite	11784
1	H5	AB-005 azepane isomer	11877
1	H6	EAM2201	11916
1	H7	UR-144 N-(5-chloropentyl) analog	11951
1	H8	THJ 018	11962
1	H9	RCS-4 M9 metabolite	11977
1	H10	RCS-4 M10 metabolite	11978
1	H11	UR-144 N-(5-bromopentyl) analog	12003
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	JWH 387	12020
2	A3	MAM2201 N-(5-chloropentyl) analog	13103
2	A4	UR-144 N-(4-chloropentyl) analog	13117
2	A5	(±)5- <i>epi</i> CP 55,940	13803
2	A6	JWH 071	13890
2	A7	A-836339	13898
2	A8	JWH 116	13977
2	A9	JWH 149	13978
2	A10	JWH 167	13979
2	A11	JWH 193	13980
2	A12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 4 of 19

Plate	Well	Contents	Item Number
2	B1	<i>Unused</i>	
2	B2	JWH 198	13981
2	B3	JWH 080	14013
2	B4	XLR11 Degradant	14055
2	B5	AM1248 azepane isomer	14068
2	B6	JWH 073 6-methoxyindole analog	14084
2	B7	BB-22	14099
2	B8	5-fluoro NNEI	14147
2	B9	AKB48 N-(4-hydroxypentyl) metabolite	14221
2	B10	5-fluoro AKB48 N-(4-hydroxypentyl) metabolite	14222
2	B11	AKB48 N-(5-hydroxypentyl) metabolite	14223
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	AKB48 N-pentanoic acid metabolite	14224
2	C3	JWH 203 N-(4-hydroxypentyl) metabolite	14227
2	C4	JWH 203 N-(5-hydroxypentyl) metabolite	14228
2	C5	JWH 203 N-pentanoic acid metabolite	14229
2	C6	UR-144 N-heptyl analog	14231
2	C7	JWH 146	14232
2	C8	MN-25	14249
2	C9	ADBICA	14293
2	C10	5-Fluoropentyl-3-pyridinoylindole (hydrochloride)	14295
2	C11	5-fluoro NNEI 2'-naphthyl isomer	14318
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	RCS-4 4-hydroxyphenyl metabolite	14345
2	D3	JWH 081 N-(4-hydroxypentyl) metabolite	14348
2	D4	JWH 081 N-pentanoic acid metabolite	14353
2	D5	XLR11 6-hydroxyindole metabolite	14354
2	D6	AM694 N-(5-hydroxypentyl) metabolite	14358
2	D7	AM694 N-pentanoic acid metabolite	14359
2	D8	BB-22 3-carboxyindole metabolite	14378
2	D9	5-fluoro PB-22 3-carboxyindole metabolite	14381
2	D10	PB-22 3-carboxyindole metabolite	14385
2	D11	PB-22 N-(4-hydroxypentyl) metabolite	14386
2	D12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 5 of 19

Plate	Well	Contents	Item Number
2	E1	<i>Unused</i>	
2	E2	PB-22 N-(5-hydroxypentyl) metabolite	14387
2	E3	PB-22 N-pentanoic acid metabolite	14389
2	E4	5-fluoro JWH 018 adamantyl analog	14432
2	E5	UR-144 N-(2-hydroxypentyl) metabolite	14470
2	E6	BB-22 3-hydroxyquinoline isomer	14505
2	E7	BB-22 4-hydroxyquinoline isomer	14506
2	E8	BB-22 5-hydroxyquinoline isomer	14507
2	E9	BB-22 6-hydroxyquinoline isomer	14508
2	E10	5-fluoro PB-22 3-hydroxyquinoline isomer	14511
2	E11	5-fluoro PB-22 4-hydroxyquinoline isomer	14512
2	E12	<i>Unused</i>	
2	F1	<i>Unused</i>	
2	F2	5-fluoro PB-22 5-hydroxyquinoline isomer	14513
2	F3	5-fluoro PB-22 6-hydroxyquinoline isomer	14514
2	F4	5-fluoro PB-22 7-hydroxyquinoline isomer	14515
2	F5	PB-22 3-hydroxyquinoline isomer	14516
2	F6	PB-22 4-hydroxyquinoline isomer	14517
2	F7	PB-22 5-hydroxyquinoline isomer	14518
2	F8	PB-22 6-hydroxyquinoline isomer	14519
2	F9	PB-22 7-hydroxyquinoline isomer	14521
2	F10	BB-22 4-hydroxyisoquinoline isomer	14541
2	F11	BB-22 5-hydroxyisoquinoline isomer	14542
2	F12	<i>Unused</i>	
2	G1	<i>Unused</i>	
2	G2	BB-22 6-hydroxyisoquinoline isomer	14543
2	G3	BB-22 7-hydroxyisoquinoline isomer	14544
2	G4	BB-22 8-hydroxyisoquinoline isomer	14545
2	G5	5-fluoro PB-22 N-(3-fluoropentyl) isomer	14547
2	G6	5-fluoro PB-22 4-hydroxyisoquinoline isomer	14549
2	G7	5-fluoro PB-22 5-hydroxyisoquinoline isomer	14550
2	G8	5-fluoro PB-22 6-hydroxyisoquinoline isomer	14551
2	G9	5-fluoro PB-22 7-hydroxyisoquinoline isomer	14552
2	G10	5-fluoro PB-22 8-hydroxyisoquinoline isomer	14553
2	G11	PB-22 5-hydroxyisoquinoline isomer	14555
2	G12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 6 of 19

Plate	Well	Contents	Item Number
2	H1	<i>Unused</i>	
2	H2	PB-22 6-hydroxyisoquinoline isomer	14556
2	H3	PB-22 8-hydroxyisoquinoline isomer	14558
2	H4	PB-22 N-(5-hydroxypentyl)-3-carboxyindole metabolite	14564
2	H5	PB-22 N-pentanoic acid-3-carboxyindole metabolite	14565
2	H6	JWH 018 N-propanoic acid metabolite	14646
2	H7	JWH 081-N-(cyclohexylmethyl) analog	14655
2	H8	NNEI 2'-naphthyl isomer	14738
2	H9	5-fluoro ADB-PINACA	14764
2	H10	5-fluoro ADBICA	14766
2	H11	1'-Naphthoyl-2-methylindole	14801
2	H12	<i>Unused</i>	
3	A1	<i>Unused</i>	
3	A2	JWH 019 N-(5-hydroxyhexyl) metabolite	14814
3	A3	5-fluoro MN-18	14816
3	A4	MN-18	14817
3	A5	THJ	14889
3	A6	5-fluoro THJ	14890
3	A7	JWH 018 8-quinolinyl carboxamide	14891
3	A8	AM2201 8-quinolinyl carboxamide	14892
3	A9	XLR12	14935
3	A10	FUB-PB-22	14949
3	A11	FDU-PB-22	14968
3	A12	<i>Unused</i>	
3	B1	<i>Unused</i>	
3	B2	NNEI	15001
3	B3	AB-PINACA N-(4-hydroxypentyl) metabolite	15049
3	B4	AB-PINACA N-(5-hydroxypentyl) metabolite	15050
3	B5	AB-PINACA N-pentanoic acid metabolite	15051
3	B6	ADB-PINACA N-(4-hydroxypentyl) metabolite	15053
3	B7	ADB-PINACA N-(5-hydroxypentyl) metabolite	15054
3	B8	ADB-PINACA pentanoic acid metabolite	15055
3	B9	ADBICA N-(5-hydroxypentyl) metabolite	15060
3	B10	ADBICA N-pentanoic acid metabolite	15061
3	B11	CI2201	15065
3	B12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 7 of 19

Plate	Well	Contents	Item Number
3	C1	<i>Unused</i>	
3	C2	JWH 018 benzimidazole analog	15074
3	C3	SDB-006	15156
3	C4	ADBICA N-(4-hydroxypentyl) metabolite	15240
3	C5	5-fluoro SDB-005	15353
3	C6	Mepirapim (hydrochloride)	15388
3	C7	SDB-005	15389
3	C8	5-fluoro SDB-006	15390
3	C9	AB-PINACA N-(2-fluoropentyl) isomer	15439
3	C10	JWH 019 N-(3-fluorohexyl) isomer	15443
3	C11	JWH 019 N-(4-fluorohexyl) isomer	15444
3	C12	<i>Unused</i>	
3	D1	<i>Unused</i>	
3	D2	AMB	15488
3	D3	AB-FUBINACA metabolite 2A	15529
3	D4	EG 018	15533
3	D5	NPB-22	15535
3	D6	5-fluoro NPB-22	15536
3	D7	PF-03550096	15620
3	D8	MMB018	15970
3	D9	MMB2201	15971
3	D10	5-chloro NNEI	16003
3	D11	M-144	16084
3	D12	<i>Unused</i>	
3	E1	<i>Unused</i>	
3	E2	MDMB-CHMINACA	16200
3	E3	PX 1	16201
3	E4	AB-CHMINACA metabolite M1B	16388
3	E5	AB-CHMINACA metabolite M2	16389
3	E6	AB-CHMINACA metabolite M3A	16391
3	E7	AB-CHMINACA metabolite M4	16393
3	E8	AB-CHMINACA metabolite M5A	16394
3	E9	MA-CHMINACA	16421
3	E10	PX 2	16434
3	E11	CBL-018	16443
3	E12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 8 of 19

Plate	Well	Contents	Item Number
3	F1	<i>Unused</i>	
3	F2	AB-CHMINACA metabolite M6	16453
3	F3	FUB-NPB-22	16883
3	F4	EG2201	16884
3	F5	Azidoindolene 1	17083
3	F6	MO-CHMINACA	17136
3	F7	5-fluoro-2-ADB-PINACA isomer 2	17169
3	F8	APP-FUBINACA	17200
3	F9	CUMYL-THPINACA	17210
3	F10	CUMYL-PICA	17211
3	F11	5-fluoro CUMYL-PICA	17212
3	F12	<i>Unused</i>	
3	G1	<i>Unused</i>	
3	G2	ADB-CHMICA	17213
3	G3	FUB-JWH 018	17332
3	G4	5-fluoro ADB-PINACA isomer 2	17396
3	G5	FDU-NNEI	17495
3	G6	JWH 307 3'-isomer	17506
3	G7	A-796260 Degradant	17515
3	G8	APP-PICA	17613
3	G9	APP-CHMINACA	17614
3	G10	AB-FUBINACA metabolite 3	17761
3	G11	MAB-CHMINACA metabolite M1	17812
3	G12	<i>Unused</i>	
3	H1	<i>Unused</i>	
3	H2	MAB-CHMINACA metabolite M3	17814
3	H3	MAB-CHMINACA metabolite M6	17815
3	H4	MAB-CHMINACA metabolite M11	17818
3	H5	5-fluoro AMB metabolite 2	17892
3	H6	5-fluoro AMB metabolite 3	17893
3	H7	AB-FUBINACA metabolite 2B	17947
3	H8	AB-CHMICA	18048
3	H9	5-chloro AKB48	18166
3	H10	3,5-AB-CHMFUPPYCA	18204
3	H11	5,3-AB-CHMFUPPYCA	18296
3	H12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 9 of 19

Plate	Well	Contents	Item Number
4	A1	<i>Unused</i>	
4	A2	5-fluoro PY-PICA	18363
4	A3	MDMB-FUBICA	18364
4	A4	ATHPINACA isomer 1	18365
4	A5	5-bromo THJ 018	18578
4	A6	5-chloro THJ 018	18581
4	A7	APINAC	18589
4	A8	5-fluoro APINAC	18592
4	A9	ADB-FUBICA	18756
4	A10	ADB-BINACA	18757
4	A11	AB-FUBICA	18758
4	A12	<i>Unused</i>	
4	B1	<i>Unused</i>	
4	B2	AB-BICA	18759
4	B3	ATHPINACA isomer 2	19040
4	B4	FUB-PB-22 3-carboxyindole metabolite	19267
4	B5	MDMB-CHMCZCA	19295
4	B6	MDMB-CHMICA metabolite M2	19316
4	B7	ADB-BICA	19729
4	B8	5-fluoro BEPIRAPIM (hydrochloride)	19730
4	B9	MDMB-FUBINACA metabolite M1	20061
4	B10	5-fluoro ADB metabolite 2	20173
4	B11	CUMYL-PICA N-pentanoic acid metabolite	20232
4	B12	<i>Unused</i>	
4	C1	<i>Unused</i>	
4	C2	MDMB-FUBICA metabolite 3	20271
4	C3	MEP-CHMICA	21652
4	C4	MEP-FUBICA	21653
4	C5	MEP-FUBINACA	21654
4	C6	CHM-122	22043
4	C7	5-fluoro MDMB-PICA metabolite 7	22095
4	C8	5-fluoro MDMB-PICA metabolite 2	22393
4	C9	7'-methoxy NABUTIE	22630
4	C10	NAPIE	22632
4	C11	4-cyano CUMYL-BUT7AICA	22818
4	C12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
4	D1	<i>Unused</i>	
4	D2	5-fluoro 7-QUPAIC	23011
4	D3	5-fluoro 7-APAICA	23021
4	D4	AB-7-FUBAICA	23059
4	D5	AB-FUBINACA 2'-indazole isomer	23065
4	D6	EADB-FUBINACA	23345
4	D7	5-fluoro- <i>tert</i> -Butylbenzyl-PINACA	23351
4	D8	5-chloro MDMB-PICA	23438
4	D9	AZEFUBIM	23653
4	D10	AEP-CHMINACA	23706
4	D11	AEP-PINACA	23707
4	D12	<i>Unused</i>	
4	E1	<i>Unused</i>	
4	E2	Ethylphenethyl-FUBICA	23764
4	E3	5-fluoro phenyl-PICA	23924
4	E4	4-cyano CUMYL-BUTCZCA	24429
4	E5	5-fluoro ethylbenzyl-PICA	24430
4	E6	Ethylbenzyl-CYBINACA	24431
4	E7	FUBIMINA N-(5-hydroxypentyl) metabolite	24468
4	E8	FUBIMINA N-pentanoic acid metabolite	24469
4	E9	MMB-FUBICA metabolite 3	24619
4	E10	ACHMINACA	25006
4	E11	MDMB-CHMCZCA metabolite M3	25064
4	E12	<i>Unused</i>	
4	F1	<i>Unused</i>	
4	F2	5-bromo APINACA	25209
4	F3	5-fluoro MDMB-7-PAICA	25286
4	F4	5-fluoro CUMYL-PeGACLONE	25435
4	F5	Isobutyl 1-pentyl-1H-indazole-3-carboxylate	25880
4	F6	MMB022	25906
4	F7	5-fluoro MPP-PICA	25916
4	F8	4-fluoro-CUMYL-5-fluoro-PICA	26131
4	F9	4-chloro CUMYL-PINACA	26132
4	F10	4-fluoro-CUMYL-5-fluoro-PINACA	26134
4	F11	4-cyano CUMYL-BUTINACA metabolite 10	26399
4	F12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
4	G1	<i>Unused</i>	
4	G2	CUMYL-CH-MeGACLONE	26432
4	G3	4-cyano ADB-BUTINACA	26506
4	G4	4-cyano MDMB-BUTINACA	26507
4	G5	5-fluoro AB-7-PAICA	26646
4	G6	APP-BUTINACA	26905
4	G7	THQ-PINACA	27081
4	G8	CUMYL-PIPETINACA (hydrochloride)	27331
4	G9	4-fluoro MDMB-BUTINACA 2'-indazole isomer	27442
4	G10	5-fluoro MDMB-PICA metabolite 8	27586
4	G11	5-fluoro MDMB-PICA metabolite 9	27587
4	G12	<i>Unused</i>	
4	H1	<i>Unused</i>	
4	H2	AB-PINACA 3-carboxyindazole metabolite	27732
4	H3	5-fluoro AB-PINACA 3-carboxyindazole metabolite	27733
4	H4	4-fluoro MDMB-BUTINACA 3-carboxyindazole metabolite	27819
4	H5	PTI-3	28238
4	H6	MDMB-3en-BUTINACA	28508
4	H7	MDMB-4en-PICA	28541
4	H8	5-fluoro MDMB-7-PAICA butanoic acid metabolite	28693
4	H9	MDMB-BUTINACA	28739
4	H10	MDMB-4en-PINACA butanoic acid metabolite	28740
4	H11	iPDMB-FUBINACA	28930
4	H12	<i>Unused</i>	
5	A1	<i>Unused</i>	
5	A2	MDMB-BUTINACA butanoic acid metabolite	28939
5	A3	4-fluoro MDMB-BUTINACA N-butanoic acid metabolite	28941
5	A4	MDMB-3en-BUTINACA butanoic acid metabolite	28942
5	A5	MDMB-4en-PICA butanoic acid metabolite	28943
5	A6	5-fluoro ADB 2'-indazole isomer	28962
5	A7	5-fluoro ADB 2'-indazole isomer butanoic acid metabolite	28963
5	A8	4-fluoro MDMB-BUTINACA 2'-isomer butanoic acid metabolite	28964
5	A9	NE-CHMIMO	29301
5	A10	MMB-4en-PINACA	29750
5	A11	5-chloro AMB-PICA	29751
5	A12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
5	B1	<i>Unused</i>	
5	B2	5-bromo AMB-PICA	29752
5	B3	CUMYL-CBMICA	30207
5	B4	ADB-BUTINACA N-butanoic acid metabolite	30294
5	B5	ADB-BUTINACA N-(4-hydroxybutyl) metabolite	30295
5	B6	MMB-4en-PINACA butanoic acid metabolite	30301
5	B7	AMB butanoic acid metabolite	30302
5	B8	CBGP	30312
5	B9	4-fluoro ABUTINACA	30724
5	B10	4-fluoro ABUTINACA N-butanoic acid metabolite	30749
5	B11	4-fluoro ABUTINACA N-(4-hydroxybutyl) metabolite	30750
5	B12	<i>Unused</i>	
5	C1	<i>Unused</i>	
5	C2	5-fluoro EMB-PICA	30769
5	C3	MMB-FUBGACLONE	30770
5	C4	4-fluoro MDMB-BUTICA butanoic acid metabolite	31106
5	C5	AFUB7AICA	31185
5	C6	MDMB-CHM7AICA	31186
5	C7	MDMB-FUB7AICA	31187
5	C8	MMB-CHM7AICA	31188
5	C9	5-fluoro MMB-P7AICA	31189
5	C10	5-fluoro MMB-PICA butanoic acid metabolite	31267
5	C11	MPP-PICA	31268
5	C12	<i>Unused</i>	
5	D1	<i>Unused</i>	
5	D2	5-fluoro MPP-PICA phenylpropionic acid metabolite	31271
5	D3	MMB-4en-PICA butanoic acid metabolite	31275
5	D4	APP-BUTINACA phenylpropanoic acid metabolite	31279
5	D5	APP-PICA phenylpropanoic acid metabolite	31280
5	D6	5-fluoro CUMYL-PeGACLONE N-(5-hydroxypentyl) metabolite	31283
5	D7	5-fluoro CUMYL-PeGACLONE N-pentanoic acid metabolite	31284
5	D8	5-fluoro CUMYL-PINACA N-pentanoic acid metabolite	31286
5	D9	4-fluoro EMB-BUTICA	31298
5	D10	Benzyl-4-cyano BUTINACA	31313
5	D11	CUMYL-CBMINACA	31325
5	D12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
5	E1	<i>Unused</i>	
5	E2	CUMYL-CB-MeGACLONE	31427
5	E3	5-fluoro EMB-PICA N-(5-hydroxypentyl) metabolite	31540
5	E4	4-fluoro MDMA-BUTICA N-(4-hydroxybutyl) metabolite	31541
5	E5	4-fluoro MDMA-BUTINACA N-(butanoic acid) 3,3-dimethylbutanoic acid metabolite	31643
5	E6	5-fluoro CUMYL-P7AICA N-pentanoic acid metabolite	31749
5	E7	4-chloro MDMA-BUTICA	31805
5	E8	CUMYL-NB-MeGACLONE	32793
5	E9	JWH 184	32794
5	E10	AFUB7AICA 7'-azaindole isomer	32989
5	E11	ADB-PHETINACA	33194
5	E12	<i>Unused</i>	
5	F1	<i>Unused</i>	
5	F2	4-cyano MMB-BUTINACA	33334
5	F3	CUMYL-NBMICA	33372
5	F4	4-fluoro MDMA-BUTINACA	33374
5	F5	MDMA-PINACA	33663
5	F6	4-cyano MMB-BUTINACA butanoic acid metabolite	33668
5	F7	AMP-4en-PINACA	33818
5	F8	ADB-HEXINACA	33820
5	F9	4-cyano MMB-BUTINACA N-butanoic acid metabolite	33835
5	F10	4-cyano MMB-BUTINACA N-butanoic acid 3-methylbutanoic acid metabolite	33836
5	F11	ADB-P7AICA	34352
5	F12	<i>Unused</i>	
5	G1	<i>Unused</i>	
5	G2	5-fluoro ADB-P7AICA	34353
5	G3	5-fluoro BZO-POXIZID	34586
5	G4	BZO-POXIZID	34795
5	G5	CUMYL-NBMINACA	34908
5	G6	BZO-CHMOXIZID	34956
5	G7	ADB-FUBIATA	35059
5	G8	MDMA-BENZICA	35135
5	G9	4-cyano AB-BUTICA	35192
5	G10	ADB-FUBIATA 3,3-dimethylbutanoic acid metabolite	35222
5	G11	MN-18 N-(5-hydroxypentyl) metabolite	35278
5	G12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
5	H1	<i>Unused</i>	
5	H2	BZO-4en-POXIZID	35503
5	H3	ADB-HEXINACA N-(6-hydroxyhexyl) metabolite	35566
5	H4	ADB-HEXINACA 3,3-dimethylbutanoic acid metabolite	35567
5	H5	ADB-HEXINACA N-hexanoic acid metabolite	35773
5	H6	ADB-BUTINAATA	36014
5	H7	2-(1-(4-Fluorobenzyl)-1H-indol-3-yl)acetic Acid	36015
5	H8	MDA-19 N-(5-hydroxyhexyl) metabolite	36039
5	H9	MDA-19 4-hydroxybenzoyl metabolite	36040
5	H10	3,5-ADB-4en-PFUPPYCA	36141
5	H11	5,3-ADB-4en-PFUPPYCA	36146
5	H12	<i>Unused</i>	
6	A1	<i>Unused</i>	
6	A2	ADB-5Br-INACA	36263
6	A3	EDMB-4en-PINACA	36444
6	A4	CH-PIATA	36558
6	A5	MDMB-5Br-INACA	36757
6	A6	ADB-5'Br-BUTINACA	36770
6	A7	CUMYL-TsINACA	36834
6	A8	MPP-BUTINACA	36852
6	A9	ADB-FUBINAATA	36853
6	A10	MMB-5Br-INACA	36883
6	A11	BZO-HEPOXIZID	36946
6	A12	<i>Unused</i>	
6	B1	<i>Unused</i>	
6	B2	AFUBIATA	36992
6	B3	ADB-IATA	37053
6	B4	CH-FUBIATA	37054
6	B5	CUMYL-INACA	37055
6	B6	CUMYL-CHSINACA	37057
6	B7	5-fluoro AMB metabolite 6	37058
6	B8	2-fluoro AMB	37319
6	B9	CHO-4'Me-5'Br-FUBOXPYRA	37468
6	B10	CH-PIATA N-pentanoic acid metabolite	37603
6	B11	ADB-5'Br-BUTINACA 3,3-dimethyl butanoic acid metabolite	37607
6	B12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
6	C1	<i>Unused</i>	
6	C2	4-fluoro AB-BUTINACA	37657
6	C3	MDMB-5'Br-BUTINACA	37698
6	C4	ADB-5'Br-DECINACA	37721
6	C5	ADB-INACA	37729
6	C6	ADB-5'Br-4en-PINACA	37741
6	C7	A-PONASA	37742
6	C8	ADB-5'Br-PINACA	37758
6	C9	5-fluoro ADB-5'Br-PINACA	37963
6	C10	JWH 181	38120
6	C11	AB-CHMIATA	38136
6	C12	<i>Unused</i>	
6	D1	<i>Unused</i>	
6	D2	MDMB-INACA 3,3-dimethylbutanoic acid metabolite	38151
6	D3	CHM-FUBIATA	38205
6	D4	CH-HEXIATA	38283
6	D5	NMDMSB	38886
6	D6	MDMB-5'Br-4en-PINACA	39488
6	D7	MDMB-5Me-INACA	39658
6	D8	MDMB-MINACA	39659
6	D9	MDMB-PICA	39729
6	D10	MDMB-PINACA	39730
6	D11	MDMB-BUTICA	40205
6	D12	<i>Unused</i>	
6	E1	<i>Unused</i>	
6	E2	HU-308	90086
6	E3	JWH 019 N-(6-hydroxyhexyl) metabolite	9000765
6	E4	JWH 250 5-hydroxyindole metabolite	9000766
6	E5	JWH 250 N-(5-hydroxypentyl) metabolite	9000767
6	E6	JWH 081 N-(5-hydroxypentyl) metabolite	9000768
6	E7	JWH 210 5-hydroxyindole metabolite	9000771
6	E8	JWH 210 N-(5-hydroxypentyl) metabolite	9000772
6	E9	JWH 018 adamantyl analog	9000799
6	E10	JWH 018 4-hydroxyindole metabolite	9000851
6	E11	JWH 018 5-hydroxyindole metabolite	9000852
6	E12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
6	F1	<i>Unused</i>	
6	F2	JWH 018 7-hydroxyindole metabolite	9000854
6	F3	JWH 018 N-(5-hydroxypentyl) metabolite	9000855
6	F4	JWH 018 N-pentanoic acid metabolite	9000856
6	F5	JWH 073 4-hydroxyindole metabolite	9000861
6	F6	JWH 073 6-hydroxyindole metabolite	9000863
6	F7	JWH 073 N-(4-hydroxybutyl) metabolite	9000865
6	F8	JWH 073 N-butanoic acid metabolite	9000866
6	F9	JWH 018 N-(2,2-dimethylpropyl) isomer	9001001
6	F10	JWH 018 N-(1,2-dimethylpropyl) isomer	9001003
6	F11	JWH 018 2'-naphthyl isomer	9001004
6	F12	<i>Unused</i>	
6	G1	<i>Unused</i>	
6	G2	JWH 018 2'-naphthyl-N-(1,2-dimethylpropyl) isomer	9001006
6	G3	JWH 018 2'-naphthyl-N-(2,2-dimethylpropyl) isomer	9001008
6	G4	JWH 073 2'-naphthyl isomer	9001014
6	G5	JWH 073 2'-naphthyl-N-(2-methylpropyl) isomer	9001015
6	G6	JWH 073 2'-naphthyl-N-(1-methylpropyl) isomer	9001016
6	G7	JWH 251 3-methylphenyl isomer	9001021
6	G8	JWH 251 4-methylphenyl isomer	9001022
6	G9	JWH 398 2-chloronaphthyl isomer	9001023
6	G10	JWH 398 5-chloronaphthyl isomer	9001025
6	G11	JWH 398 6-chloronaphthyl isomer	9001026
6	G12	<i>Unused</i>	
6	H1	<i>Unused</i>	
6	H2	JWH 398 7-chloronaphthyl isomer	9001027
6	H3	AM2201 N-(3-fluoropentyl) isomer	9001030
6	H4	JWH 122 3-methylnaphthyl isomer	9001033
6	H5	JWH 122 6-methylnaphthyl isomer	9001035
6	H6	JWH 122 7-methylnaphthyl isomer	9001036
6	H7	JWH 122 8-methylnaphthyl isomer	9001037
6	H8	JWH 210 2-ethylnaphthyl isomer	9001038
6	H9	JWH 210 8-ethylnaphthyl isomer	9001043
6	H10	JWH 081 2-methoxynaphthyl isomer	9001044
6	H11	JWH 081 3-methoxynaphthyl isomer	9001045
6	H12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
7	A1	<i>Unused</i>	
7	A2	JWH 081 6-methoxynaphthyl isomer	9001047
7	A3	JWH 081 8-methoxynaphthyl isomer	9001049
7	A4	AM1220	9001055
7	A5	JWH 022	9001056
7	A6	JWH 073 4-methylnaphthyl analog	9001076
7	A7	JWH 018 adamantyl carboxamide	9001193
7	A8	JWH 072	9001201
7	A9	JWH 424	9001203
7	A10	JWH 180	9001205
7	A11	MAM2201	9001219
7	A12	<i>Unused</i>	
7	B1	<i>Unused</i>	
7	B2	UR-144 N-(5-methylhexyl) analog	9001439
7	B3	UR-144 Degradant N-pentanoic acid metabolite	9001453
7	B4	MN-25-2-methyl derivative	9001455
7	B5	AB-FUBINACA 3-fluorobenzyl isomer	9001523
7	B6	AB-FUBINACA 2-fluorobenzyl isomer	9001524
7	B7	ADB-PINACA isomer 1	9001525
7	B8	ADB-PINACA isomer 2	9001526
7	B9	ADB-PINACA isomer 3	9001527
7	B10	AB-FUBINACA isomer 1	9001528
7	B11	AB-FUBINACA isomer 2	9001529
7	B12	<i>Unused</i>	
7	C1	<i>Unused</i>	
7	C2	AB-FUBINACA isomer 5	9001546
7	C3	5-fluoro ABICA	9001555
7	C4	F2201	9001580
7	C5	5-chloro AB-PINACA	9001857
7	C6	PTI-1 (hydrochloride)	9001948
7	C7	(1H-Indol-3-yl)(2,2,3,3-tetramethylcyclopropyl) methanone	9001966
7	C8	ADB-PINACA isomer 4	9001994
7	C9	4-fluoro ADB	9002717
7	C10	3-fluoro AMB	9002737
7	C11	4-fluoro AMB	9002738
7	C12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 18 of 19

Plate	Well	Contents	Item Number
7	D1	Unused	
7	D2	2-fluoro NNEI	9002739
7	D3	3-fluoro NNEI	9002740
7	D4	Ethylbenzyl-PeGACLONE	9003353
7	D5	THJ2201 N-(5-hydroxypentyl) metabolite	9003736
7	D6	JWH 133	10005428
7	D7	HU-211	10006350
7	D8	Pravadoline	10006973
7	D9	JWH 015	10009018
7	D10	STS-135	ISO00058
7	D11	Unused	
7	D12	Unused	
7	E1	Unused	
7	E2	Unused	
7	E3	Unused	
7	E4	Unused	
7	E5	Unused	
7	E6	Unused	
7	E7	Unused	
7	E8	Unused	
7	E9	Unused	
7	E10	Unused	
7	E11	Unused	
7	E12	Unused	
7	F1	Unused	
7	F2	Unused	
7	F3	Unused	
7	F4	Unused	
7	F5	Unused	
7	F6	Unused	
7	F7	Unused	
7	F8	Unused	
7	F9	Unused	
7	F10	Unused	
7	F11	Unused	
7	F12	Unused	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

[illegible]

WWW.CAYMANCHEM.COM