

PRODUCT INFORMATION



Autophagy-focused Screening Library (96-Well)

Item No. 23537 • Batch No. 0517796

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

Page 1 of 6

| Plate | Well | Contents | Item Number |
|-------|------|-------------------------|-------------|
| 1 | A1 | <i>Unused</i> | |
| 1 | A2 | SB939 | 10443 |
| 1 | A3 | PCI 34051 | 10444 |
| 1 | A4 | Sirtinol | 10523 |
| 1 | A5 | Scriptaid | 10572 |
| 1 | A6 | IU1 | 10617 |
| 1 | A7 | Pifithrin- μ | 10748 |
| 1 | A8 | Piperlongumine | 11006 |
| 1 | A9 | Delphinidin (chloride) | 11012 |
| 1 | A10 | Forskolin | 11018 |
| 1 | A11 | SMI-4a | 11029 |
| 1 | A12 | <i>Unused</i> | |
| 1 | B1 | <i>Unused</i> | |
| 1 | B2 | ITF 2357 | 11045 |
| 1 | B3 | PFI-1 | 11155 |
| 1 | B4 | Sodium 4-Phenylbutyrate | 11323 |
| 1 | B5 | YM-155 | 11490 |
| 1 | B6 | Obatoclox (mesylate) | 11499 |
| 1 | B7 | ABT-737 | 11501 |
| 1 | B8 | ABT-888 (hydrochloride) | 11505 |
| 1 | B9 | IOX2 | 11573 |
| 1 | B10 | Temsirolimus | 11590 |
| 1 | B11 | Everolimus | 11597 |
| 1 | B12 | <i>Unused</i> | |
| 1 | C1 | <i>Unused</i> | |
| 1 | C2 | AZD 1152-HQPA | 11602 |
| 1 | C3 | Tamoxifen (citrate) | 11629 |
| 1 | C4 | Gemcitabine | 11690 |
| 1 | C5 | Vinblastine (sulfate) | 11762 |
| 1 | C6 | Vincristine (sulfate) | 11764 |
| 1 | C7 | Dioscin | 11834 |
| 1 | C8 | Brefeldin A | 11861 |
| 1 | C9 | A-769662 | 11900 |
| 1 | C10 | Dorsomorphin | 11967 |
| 1 | C11 | Rottlerin | 12006 |
| 1 | C12 | <i>Unused</i> | |

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, 05/08/2026

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 6

| Plate | Well | Contents | Item Number |
|-------|------|-------------------------------------------|-------------|
| 1 | D1 | <i>Unused</i> | |
| 1 | D2 | CI-994 | 12084 |
| 1 | D3 | 2-Methoxyestradiol | 13021 |
| 1 | D4 | Valproic Acid (sodium salt) | 13033 |
| 1 | D5 | STF-62247 | 13084 |
| 1 | D6 | Tenovin-1 | 13085 |
| 1 | D7 | Metformin (hydrochloride) | 13118 |
| 1 | D8 | M 344 | 13174 |
| 1 | D9 | 3-Methyladenine | 13242 |
| 1 | D10 | Tamoxifen | 13258 |
| 1 | D11 | MS-275 | 13284 |
| 1 | D12 | <i>Unused</i> | |
| 1 | E1 | <i>Unused</i> | |
| 1 | E2 | FK-866 | 13287 |
| 1 | E3 | Pifithrin- α | 13326 |
| 1 | E4 | Rapamycin | 13346 |
| 1 | E5 | Geldanamycin | 13355 |
| 1 | E6 | MK-0457 | 13600 |
| 1 | E7 | ZM 447439 | 13601 |
| 1 | E8 | MLN8237 | 13602 |
| 1 | E9 | WYE-354 | 13604 |
| 1 | E10 | PP242 | 13643 |
| 1 | E11 | Nocodazole | 13857 |
| 1 | E12 | <i>Unused</i> | |
| 1 | F1 | <i>Unused</i> | |
| 1 | F2 | JNJ-26481585 (hydrochloride) | 14088 |
| 1 | F3 | Temozolomide | 14163 |
| 1 | F4 | Torin 2 | 14185 |
| 1 | F5 | Chloroquine (phosphate) | 14194 |
| 1 | F6 | 2-deoxy-D-Glucose | 14325 |
| 1 | F7 | Acridine Orange | 14338 |
| 1 | F8 | Nimodipine | 14573 |
| 1 | F9 | GSK2578215A | 14603 |
| 1 | F10 | (\pm)-Thalidomide | 14610 |
| 1 | F11 | Cyclic Pifithrin- α (hydrobromide) | 14748 |
| 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 6

| Plate | Well | Contents | Item Number |
|-------|------|-------------------------------------|-------------|
| 1 | G1 | <i>Unused</i> | |
| 1 | G2 | Nafamostat (mesylate) | 14837 |
| 1 | G3 | Amlodipine | 14838 |
| 1 | G4 | Loperamide (hydrochloride) | 14875 |
| 1 | G5 | Omeprazole | 14880 |
| 1 | G6 | SNS-314 (mesylate) | 14974 |
| 1 | G7 | Azithromycin | 15004 |
| 1 | G8 | Doxorubicin (hydrochloride) | 15007 |
| 1 | G9 | PF-04691502 | 15017 |
| 1 | G10 | Trifluoperazine (hydrochloride) | 15068 |
| 1 | G11 | Amiodarone (hydrochloride) | 15213 |
| 1 | G12 | <i>Unused</i> | |
| 1 | H1 | <i>Unused</i> | |
| 1 | H2 | P005091 | 15224 |
| 1 | H3 | WP1130 | 15227 |
| 1 | H4 | Roxadustat | 15294 |
| 1 | H5 | DBeQ | 15318 |
| 1 | H6 | Taurolidine | 15374 |
| 1 | H7 | LDN-57444 | 15396 |
| 1 | H8 | Carbamazepine | 15422 |
| 1 | H9 | Ranolazine (hydrochloride) | 15604 |
| 1 | H10 | <i>Unused</i> | |
| 1 | H11 | Clonidine (hydrochloride) | 15949 |
| 1 | H12 | <i>Unused</i> | |
| 2 | A1 | <i>Unused</i> | |
| 2 | A2 | <i>p</i> -nitro-Pifithrin- α | 16209 |
| 2 | A3 | GNE-0877 | 16251 |
| 2 | A4 | GNE-7915 | 16252 |
| 2 | A5 | JNJ-26854165 | 16259 |
| 2 | A6 | MC 1568 | 16265 |
| 2 | A7 | ONX 0912 | 16269 |
| 2 | A8 | ONX 0914 | 16271 |
| 2 | A9 | XL147 analog 2 | 16299 |
| 2 | A10 | SGI-1776 | 16423 |
| 2 | A11 | CUDC-101 | 16426 |
| 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 6

| Plate | Well | Contents | Item Number |
|-------|------|----------------------------------------|-------------|
| 2 | B1 | <i>Unused</i> | |
| 2 | B2 | LAQ824 | 16427 |
| 2 | B3 | Bupivacaine | 16618 |
| 2 | B4 | AZD 8055 | 16978 |
| 2 | B5 | SAR405 | 16979 |
| 2 | B6 | Perhexiline (maleate) | 16982 |
| 2 | B7 | PIK-III | 17002 |
| 2 | B8 | VLX600 | 17192 |
| 2 | B9 | DL- <i>threo</i> -PPMP (hydrochloride) | 17236 |
| 2 | B10 | Verteporfin | 17334 |
| 2 | B11 | ML-240 | 17373 |
| 2 | B12 | <i>Unused</i> | |
| 2 | C1 | <i>Unused</i> | |
| 2 | C2 | OSI-027 | 17379 |
| 2 | C3 | VPS34-IN-1 | 17392 |
| 2 | C4 | (±)-Felodipine | 17535 |
| 2 | C5 | Isradipine | 17536 |
| 2 | C6 | Nitrendipine | 17549 |
| 2 | C7 | Carfilzomib | 17554 |
| 2 | C8 | NMS-873 | 17674 |
| 2 | C9 | Quinacrine analog 34 | 17733 |
| 2 | C10 | CE3F4 | 17767 |
| 2 | C11 | SMER28 | 17768 |
| 2 | C12 | <i>Unused</i> | |
| 2 | D1 | <i>Unused</i> | |
| 2 | D2 | Spatin-1 | 17769 |
| 2 | D3 | Hydroxychloroquine (sulfate) | 17911 |
| 2 | D4 | MHY1485 | 18008 |
| 2 | D5 | LRRK2-IN-1 | 18094 |
| 2 | D6 | CCT128930 | 18194 |
| 2 | D7 | Mocetinostat | 18287 |
| 2 | D8 | PF-06447475 | 18375 |
| 2 | D9 | MLN9708 | 18386 |
| 2 | D10 | Danusertib | 18387 |
| 2 | D11 | SBI-0206965 | 18477 |
| 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 6

| Plate | Well | Contents | Item Number |
|-------|------|------------------------------|-------------|
| 2 | E1 | <i>Unused</i> | |
| 2 | E2 | GSK621 | 19155 |
| 2 | E3 | MK-5108 | 19167 |
| 2 | E4 | CPTH6 (hydrobromide) | 19828 |
| 2 | E5 | Chrysophanol | 19870 |
| 2 | E6 | MRT68921 | 19905 |
| 2 | E7 | PCI 24781 | 20059 |
| 2 | E8 | Sulfacetamide (sodium salt) | 20377 |
| 2 | E9 | Flunarizine (hydrochloride) | 20385 |
| 2 | E10 | L-Tetrahydropalmatine | 20535 |
| 2 | E11 | PP121 | 21196 |
| 2 | E12 | <i>Unused</i> | |
| 2 | F1 | <i>Unused</i> | |
| 2 | F2 | CYC-116 | 21205 |
| 2 | F3 | Dorsomorphin (hydrochloride) | 21207 |
| 2 | F4 | Oleuropein | 21220 |
| 2 | F5 | Nilvadipine | 21243 |
| 2 | F6 | Azelnidipine | 21249 |
| 2 | F7 | ACY-1215 | 21531 |
| 2 | F8 | Aurora A Inhibitor I | 21600 |
| 2 | F9 | ZLN024 | 21644 |
| 2 | F10 | Ro 3280 | 21669 |
| 2 | F11 | NVP-BGT226 | 22142 |
| 2 | F12 | <i>Unused</i> | |
| 2 | G1 | <i>Unused</i> | |
| 2 | G2 | Divalproex (sodium salt) | 22960 |
| 2 | G3 | Clevidipine Butyrate | 23025 |
| 2 | G4 | GNE-9605 | 23446 |
| 2 | G5 | Parthenolide | 70080 |
| 2 | G6 | Aspirin | 70260 |
| 2 | G7 | <i>trans</i> -Resveratrol | 70675 |
| 2 | G8 | LY294002 | 70920 |
| 2 | G9 | Trichostatin A | 89730 |
| 2 | G10 | (E/Z)-Belinostat | 10004279 |
| 2 | G11 | Y-27632 (hydrochloride) | 10005583 |
| 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 6

| Plate | Well | Contents | Item Number |
|-------|------|--------------------------|-------------|
| 2 | H1 | <i>Unused</i> | |
| 2 | H2 | RITA | 10006426 |
| 2 | H3 | PI-103 | 10009209 |
| 2 | H4 | EX-527 | 10009798 |
| 2 | H5 | SAHA | 10009929 |
| 2 | H6 | BAY 11-7082 | 10010266 |
| 2 | H7 | SP600125 | 10010466 |
| 2 | H8 | HA-1077 (hydrochloride) | 10010559 |
| 2 | H9 | Wortmannin | 10010591 |
| 2 | H10 | SRT 1720 (hydrochloride) | 10011020 |
| 2 | H11 | (S)-MG132 | 10012628 |
| 2 | 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