

PRODUCT INFORMATION



Prostaglandin Screening Library I (96-Well)

Item No. 10501 • Batch No. 0609348

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 | Unused | |
| 1 | A2 | Prostaglandin F _{2α} -1-glyceryl ester | 10139 |
| 1 | A3 | Prostaglandin F _{2α} serinol amide | 10194 |
| 1 | A4 | Tafluprost ethyl ester | 11612 |
| 1 | A5 | 13,14-dihydro-16,16-difluoro Prostaglandin F _{2α} | 13609 |
| 1 | A6 | 13,14-epoxy Fluprostenol isopropyl ester | 13679 |
| 1 | A7 | Prostaglandin F _{1α} | 15010 |
| 1 | A8 | Prostaglandin F _{1α} Alcohol | 15020 |
| 1 | A9 | 2,3-dinor-6-keto Prostaglandin F _{1α} (sodium salt) | 15120 |
| 1 | A10 | 3-methoxy Prostaglandin F _{1α} | 15170 |
| 1 | A11 | 6-keto Prostaglandin F _{1α} | 15210 |
| 1 | A12 | Unused | |
| 1 | B1 | Unused | |
| 1 | B2 | 6,15-diketo-13,14-dihydro Prostaglandin F _{1α} | 15270 |
| 1 | B3 | 8-iso Prostaglandin F _{1α} | 15350 |
| 1 | B4 | 8-iso Prostaglandin F _{1β} | 15370 |
| 1 | B5 | Prostaglandin F _{1β} | 15410 |
| 1 | B6 | 11β-Prostaglandin F _{1β} | 15450 |
| 1 | B7 | 11-deoxy Prostaglandin F _{1β} | 15510 |
| 1 | B8 | 11-deoxy Prostaglandin F _{1α} | 15512 |
| 1 | B9 | 13,14-dihydro Prostaglandin F _{1α} | 15610 |
| 1 | B10 | 13,14-dihydro-15-keto Prostaglandin F _{1α} | 15670 |
| 1 | B11 | 15-keto Prostaglandin F _{1α} | 15710 |
| 1 | B12 | Unused | |
| 1 | C1 | Unused | |
| 1 | C2 | 17,20-dimethyl Prostaglandin F _{1α} | 15790 |
| 1 | C3 | 17-trifluoromethylphenyl-13,14-dihydro trinor Prostaglandin F _{1α} | 15895 |
| 1 | C4 | Prostaglandin F _{2α} | 16010 |
| 1 | C5 | Prostaglandin F _{2α} methyl ester | 16011 |
| 1 | C6 | Prostaglandin F _{2α} Alcohol | 16012 |
| 1 | C7 | Prostaglandin F _{2α} Ethanolamide | 16013 |
| 1 | C8 | Prostaglandin F _{2α} Alcohol methyl ether | 16014 |
| 1 | C9 | Prostaglandin F _{2α} ethyl amide | 16016 |
| 1 | C10 | Prostaglandin F _{2α} diethyl amide | 16023 |
| 1 | C11 | Prostaglandin F _{2α} isopropyl ester | 16030 |
| 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, 04/21/2021

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 | Prostaglandin F _{2α} dimethyl amide | 16032 |
| 1 | D3 | Prostaglandin F _{2α} dimethyl amine | 16033 |
| 1 | D4 | 5- <i>trans</i> Prostaglandin F _{2α} | 16210 |
| 1 | D5 | 8- <i>iso</i> Prostaglandin F _{2α} | 16350 |
| 1 | D6 | 8- <i>iso</i> Prostaglandin F _{2β} | 16370 |
| 1 | D7 | 8- <i>iso</i> -15(R)-Prostaglandin F _{2α} | 16395 |
| 1 | D8 | Prostaglandin F _{2β} | 16410 |
| 1 | D9 | 16,16-dimethyl Prostaglandin F _{2β} | 16470 |
| 1 | D10 | 11-deoxy Prostaglandin F _{2α} | 16500 |
| 1 | D11 | 11-deoxy Prostaglandin F _{2β} | 16510 |
| 1 | D12 | <i>Unused</i> | |
| 1 | E1 | <i>Unused</i> | |
| 1 | E2 | 11β-Prostaglandin F _{2α} | 16520 |
| 1 | E3 | 11β-Prostaglandin F _{2α} Ethanolamide | 16522 |
| 1 | E4 | 13,14-dihydro Prostaglandin F _{2α} | 16660 |
| 1 | E5 | 13,14-dihydro-15-keto Prostaglandin F _{2α} | 16670 |
| 1 | E6 | 15-keto Prostaglandin F _{2α} | 16720 |
| 1 | E7 | 15(R)-15-methyl Prostaglandin F _{2α} | 16730 |
| 1 | E8 | 15(R)-15-methyl Prostaglandin F _{2α} methyl ester | 16734 |
| 1 | E9 | 15(R)-Prostaglandin F _{2α} | 16740 |
| 1 | E10 | 15(S)-15-methyl Prostaglandin F _{2α} | 16743 |
| 1 | E11 | 15(S)-15-methyl Prostaglandin F _{2α} methyl ester | 16744 |
| 1 | E12 | <i>Unused</i> | |
| 1 | F1 | <i>Unused</i> | |
| 1 | F2 | 16,16-dimethyl Prostaglandin F _{2α} | 16750 |
| 1 | F3 | 16-phenoxy tetranor Prostaglandin F _{2α} | 16760 |
| 1 | F4 | (+)-Cloprostenol (sodium salt) | 16766 |
| 1 | F5 | Fluprostenol | 16768 |
| 1 | F6 | Fluprostenol isopropyl ester | 16769 |
| 1 | F7 | 5- <i>trans</i> Fluprostenol | 16776 |
| 1 | F8 | 5- <i>trans</i> Fluprostenol isopropyl ester | 16777 |
| 1 | F9 | 15-keto Fluprostenol | 16785 |
| 1 | F10 | 15-keto Fluprostenol isopropyl ester | 16786 |
| 1 | F11 | 15(S)-Fluprostenol | 16787 |
| 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 | 15(S)-Fluprostenol isopropyl ester | 16788 |
| 1 | G3 | 17-phenyl trinor Prostaglandin F _{2α} | 16810 |
| 1 | G4 | Latanoprost (free acid) | 16811 |
| 1 | G5 | Latanoprost | 16812 |
| 1 | G6 | 5- <i>trans</i> Latanoprost | 16813 |
| 1 | G7 | 15(R)-17-phenyl trinor Prostaglandin F _{2α} | 16814 |
| 1 | G8 | 15-keto Latanoprost (free acid) | 16815 |
| 1 | G9 | 15-keto Latanoprost | 16816 |
| 1 | G10 | 15-keto-17-phenyl trinor Prostaglandin F _{2α} | 16817 |
| 1 | G11 | 15(S)-Latanoprost | 16818 |
| 1 | G12 | <i>Unused</i> | |
| 1 | H1 | <i>Unused</i> | |
| 1 | H2 | 17-phenyl trinor Prostaglandin F _{2α} ethyl amide | 16820 |
| 1 | H3 | 17-phenyl trinor Prostaglandin F _{2α} amide | 16821 |
| 1 | H4 | Latanoprost ethyl amide | 16822 |
| 1 | H5 | 17-phenyl trinor Prostaglandin F _{2α} diethyl amide | 16823 |
| 1 | H6 | 17-phenyl trinor Prostaglandin F _{2α} isopropyl ester | 16824 |
| 1 | H7 | 15(R)-17-phenyl trinor Prostaglandin F _{2α} isopropyl ester | 16825 |
| 1 | H8 | 17-trifluoromethylphenyl trinor Prostaglandin F _{2α} | 16890 |
| 1 | H9 | 17-trifluoromethylphenyl-13,14-dihydro trinor Prostaglandin F _{2α} | 16895 |
| 1 | H10 | 20-ethyl Prostaglandin F _{2α} | 16940 |
| 1 | H11 | 15-keto Cloprostenol isopropyl ester | 24929 |
| 1 | H12 | <i>Unused</i> | |
| 2 | A1 | <i>Unused</i> | |
| 2 | A2 | (+)-5- <i>trans</i> Cloprostenol isopropyl ester | 24930 |
| 2 | A3 | 17-phenyl trinor Prostaglandin F _{2α} dimethyl amide | 25241 |
| 2 | A4 | 17-phenyl trinor Prostaglandin F _{2α} isopropyl amide | 25242 |
| 2 | A5 | Latanoprost amide | 25245 |
| 2 | A6 | Latanoprost dimethyl amide | 25247 |
| 2 | A7 | 1a-monohomo-17-phenyl trinor Prostaglandin F _{2α} | 25392 |
| 2 | A8 | 17-phenyl trinor Prostaglandin F _{2α} glycnamide | 25470 |
| 2 | A9 | 17-phenyl trinor Prostaglandin F _{2α} glycnamide methyl ester | 25471 |
| 2 | A10 | 15(R)-Prostaglandin F _{1α} | 29549 |
| 2 | A11 | 17-phenyl trinor Prostaglandin F _{2α} cyclohexyl amide | 9000686 |
| 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 | Tafluprost ethyl amide | 9000843 |
| 2 | B3 | Fluprostenol serinol amide | 10004236 |
| 2 | B4 | 17-phenyl trinor Prostaglandin F _{2α} serinol amide | 10004237 |
| 2 | B5 | (+)-5- <i>trans</i> Cloprostenol | 10004970 |
| 2 | B6 | Tafluprost (free acid) | 10005439 |
| 2 | B7 | Tafluprost | 10005440 |
| 2 | B8 | 8-iso Prostaglandin F _{2α} Ethanolamide | 10005764 |
| 2 | B9 | <i>ent</i> -Prostaglandin F _{2α} | 10008122 |
| 2 | B10 | 15(R)-17-phenyl trinor Prostaglandin F _{2α} ethyl amide | 10008127 |
| 2 | B11 | 5- <i>trans</i> -17-phenyl trinor Prostaglandin F _{2α} ethyl amide | 10008132 |
| 2 | B12 | <i>Unused</i> | |
| 2 | C1 | <i>Unused</i> | |
| 2 | C2 | 8- <i>iso</i> -17-phenyl trinor Prostaglandin F _{2α} | 10008435 |
| 2 | C3 | 8- <i>iso</i> -17-phenyl trinor Prostaglandin F _{2β} | 10008436 |
| 2 | C4 | 8- <i>iso</i> -15-keto Prostaglandin F _{2β} | 10008539 |
| 2 | C5 | 16-phenoxy Prostaglandin F _{2α} ethyl amide | 10009875 |
| 2 | C6 | (+)-Cloprostenol isopropyl ester | 10010016 |
| 2 | C7 | 17-trifluoromethylphenyl trinor Prostaglandin F _{2α} ethyl amide | 10010061 |
| 2 | C8 | 17-trifluoromethylphenyl trinor Prostaglandin F _{2α} isopropyl ester | 10010062 |
| 2 | C9 | 16-phenoxy tetranor Prostaglandin F _{2α} methyl ester | 10010102 |
| 2 | C10 | 16-phenoxy tetranor Prostaglandin F _{2α} isopropyl ester | 10010103 |
| 2 | C11 | 17-phenyl trinor Prostaglandin F _{2α} methyl ester | 10010110 |
| 2 | C12 | <i>Unused</i> | |
| 2 | D1 | <i>Unused</i> | |
| 2 | D2 | 17-trifluoromethylphenyl trinor Prostaglandin F _{2α} methyl ester | 10010111 |
| 2 | D3 | (+)-Cloprostenol methyl ester | 10010115 |
| 2 | D4 | Fluprostenol methyl ester | 10010151 |
| 2 | D5 | 17-phenyl trinor Prostaglandin F _{2α} methyl amide | 10010351 |
| 2 | D6 | 15-keto-17-phenyl trinor Prostaglandin F _{2α} ethyl amide | 10010405 |
| 2 | D7 | Fluprostenol methyl amide | 10010406 |
| 2 | D8 | (+)-Cloprostenol methyl amide | 10010495 |
| 2 | D9 | 16-phenoxy tetranor Prostaglandin F _{2α} methyl amide | 10010562 |
| 2 | D10 | 17-phenyl trinor Prostaglandin F _{2α} cyclopropyl amide | 10010605 |
| 2 | D11 | 17-phenoxy trinor Prostaglandin F _{2α} isopropyl ester | 10010723 |
| 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 | 17-phenoxy trinor Prostaglandin F _{2α} ethyl amide | 10010742 |
| 2 | E3 | 17-phenyl trinor Prostaglandin F _{2α} cyclopropyl methyl amide | 10010810 |
| 2 | E4 | 17-phenoxy trinor Prostaglandin F _{2α} | 10010839 |
| 2 | E5 | <i>Unused</i> | |
| 2 | E6 | <i>Unused</i> | |
| 2 | E7 | <i>Unused</i> | |
| 2 | E8 | <i>Unused</i> | |
| 2 | E9 | <i>Unused</i> | |
| 2 | E10 | <i>Unused</i> | |
| 2 | E11 | <i>Unused</i> | |
| 2 | E12 | <i>Unused</i> | |
| 2 | F1 | <i>Unused</i> | |
| 2 | F2 | <i>Unused</i> | |
| 2 | F3 | <i>Unused</i> | |
| 2 | F4 | <i>Unused</i> | |
| 2 | F5 | <i>Unused</i> | |
| 2 | F6 | <i>Unused</i> | |
| 2 | F7 | <i>Unused</i> | |
| 2 | F8 | <i>Unused</i> | |
| 2 | F9 | <i>Unused</i> | |
| 2 | F10 | <i>Unused</i> | |
| 2 | F11 | <i>Unused</i> | |
| 2 | F12 | <i>Unused</i> | |
| 2 | G1 | <i>Unused</i> | |
| 2 | G2 | <i>Unused</i> | |
| 2 | G3 | <i>Unused</i> | |
| 2 | G4 | <i>Unused</i> | |
| 2 | G5 | <i>Unused</i> | |
| 2 | G6 | <i>Unused</i> | |
| 2 | G7 | <i>Unused</i> | |
| 2 | G8 | <i>Unused</i> | |
| 2 | G9 | <i>Unused</i> | |
| 2 | G10 | <i>Unused</i> | |
| 2 | G11 | <i>Unused</i> | |
| 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 | Unused | |
| 2 | H2 | Unused | |
| 2 | H3 | Unused | |
| 2 | H4 | Unused | |
| 2 | H5 | Unused | |
| 2 | H6 | Unused | |
| 2 | H7 | Unused | |
| 2 | H8 | Unused | |
| 2 | H9 | Unused | |
| 2 | H10 | Unused | |
| 2 | H11 | Unused | |
| 2 | H12 | Unused | |

*Primary Target is based on available IC_{50} and/or K_i values cited in the literature

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