

Stauprimide

Primes stem cells for differentiation

Stauprimide dramatically primes embryonic stem cells for differentiation. Its molecular target is NME2, a c-Myc-activating transcription factor. Stauprimide inhibits nuclear localization of NME2 thereby downregulating c-Myc.

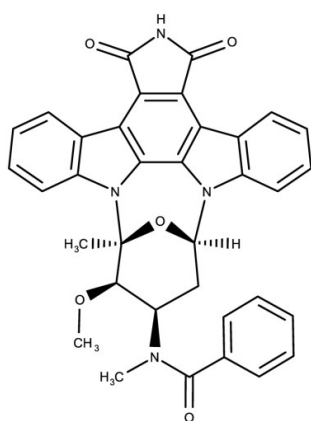
Ordering Information

[Order Online »](#)

ENZ-CHM466-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-Benzoyl-7-oxostaurosporine

Appearance Orange solid.

CAS 154589-96-5

Formula $C_{35}H_{28}N_4O_5$

Identity Determined by NMR.

MW 584.64

Purity ≥98% (TLC)

Solubility Soluble in DMSO (20 mg/ml).

Last modified: February 27, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)