

eFT508

Potent MNK1/2 inhibitor

eFT508 is a potent and highly selective inhibitor of the mitogen-activated protein kinase interacting kinases 1 and 2 (MNK1/2), key regulators of mRNA translation protein eukaryotic initiation factor eIF4E.1 It potently inhibits phosphorylation of eIF4E in multiple solid and hematological tumor lines leading to blockade of pro-inflammatory and pro-tumorigenic cytokine production. Displays inhibitory effects in various cancers models including acute myeloid leukemia, gastric, and breast. eFT-508 has also shown efficacy in treating neuropathic pain via the MNK-eIF4E pathway.

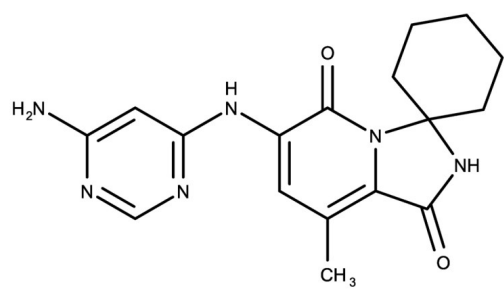
Ordering Information

[Order Online »](#)

ENZ-CHM448-0025	25mg
-----------------	------

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 6-[(6-Aminopyrimidin-4-yl)amino]-8-methylspiro[2H-imidazo[1,5-a]pyridine-3,1'-cyclohexane]-1,5-dione

Appearance Pale yellow solid.

CAS 1849590-01-7

Couple Type Inhibitor

Formula $C_{17}H_{20}N_6O_2$

Identity Determined by NMR.

MW 340.39

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (2 mg/ml with warming).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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