

PRODUCT INFORMATION

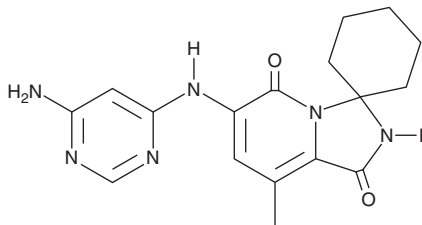


eFT508

Item No. 21957

CAS Registry No.: 1849590-01-7
Formal Name: 6'-[(6-amino-4-pyrimidinyl)amino]-8'-methyl-spiro[cyclohexane-1,3'(2'H)-imidazo[1,5-a]pyridine]-1',5'-dione

MF: C₁₇H₂₀N₆O₂
FW: 340.4
Purity: ≥98%
UV/Vis.: λ_{max}: 237, 357 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

eFT508 is supplied as a crystalline solid. A stock solution may be made by dissolving the eFT508 in the solvent of choice, which should be purged with an inert gas. eFT508 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of eFT508 in these solvents is approximately 1 and 0.5 mg/ml, respectively.

Description

eFT508 is a potent inhibitor of MAPK interacting protein kinases (MNK) 1 and 2 (IC₅₀s = <10 nM).¹ It inhibits phosphorylation of eukaryotic initiation factor 4E (eIF4E) with an IC₅₀ value of <50 nM in a cellular signaling assay. eFT508 has a half-life of 5.3 hours in rats.

Reference

1. Reich, S.H., Sprengeler, P.A., Webber, S.E., *et al.* MNK inhibitors and methods related thereto. *US 2015/0376181 A1* (2015).

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, 12/01/2022

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