

Trametinib

Potent MEK1/2 inhibitor

Trametinib is a highly specific and potent MEK1/2 inhibitor with IC_{50} of 0.92 nM/1.8 nM in cell-free assays.

Citations: 1

[View Online »](#)

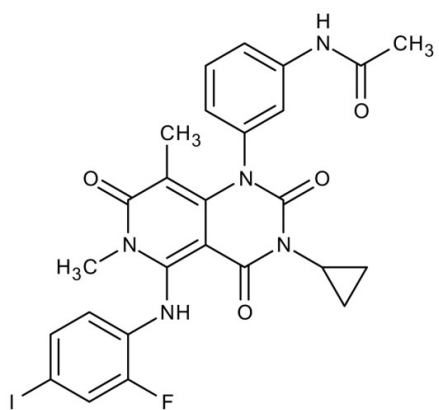
Ordering Information

[Order Online »](#)

ENZ-CHM239-0010	10mg
-----------------	------

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance White solid.

CAS 871700-17-3

Couple Target MEK

Couple Type Inhibitor

Formula $C_{26}H_{23}FIN_5O_4$

MW 615.4

Purity ≥99% (HPLC)

Solubility Soluble in DMSO (20 mg/mL).

Last modified: May 29, 2024



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-
be@enzolifesciences.com](mailto:info-be@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)