

JK 184

Hedgehog (Hh) pathway blocker.

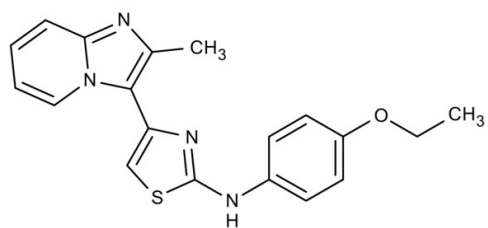
Ordering Information

[Order Online »](#)

ENZ-CHM129-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(4-ethoxyphenyl)-4-(2-methylimidazo[1,2-a]pyridin-3-yl)-1,3-thiazol-2-amine, Hh signaling antagonist VII
Appearance	Pink powder solid.
CAS	315703-52-7
Couple Target	Alcohol dehydrogenase, Sonic hedgehog
Couple Type	Inhibitor
Formula	$C_{19}H_{18}N_4OS$
MW	350.4
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO or methanol.

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-
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)