

HET-0016

Cytochrome P450 inhibitor

HET-0016 is a potent and selective inhibitor of the biosynthesis of 20-HETE acting via inhibition of CYP4A. It inhibits cerebrovascular constriction induced by Ca^{2+} or arachidonate and inhibits angiogenesis and endothelial cell sprouting in various models.

Citations: 5

[View Online »](#)

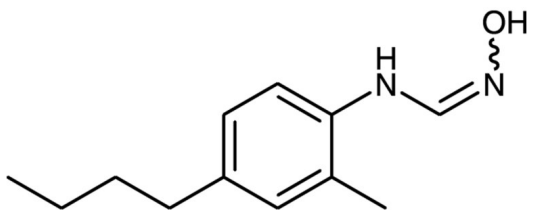
Ordering Information

[Order Online »](#)

BML-EI350-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	N-Hydroxy-N'-(4-butyl-2-methylphenyl)formamidine
Appearance	Light pink solid.
CAS	339068-25-6
Couple Target	Cytochrome P450
Couple Type	Inhibitor
Formula	$C_{12}H_{18}N_2O$
MW	206.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (6 mg/ml)

Last modified: May 29, 2024



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