

SB 590885 . dihydrochloride

Potent B-Raf inhibitor.

Selective for B-Raf against 46 other kinases. SB 590885 dihydrochloride decreases anchorage-independent growth of melanoma cell lines. Inhibits ERK phosphorylation and proliferation in tumor cells expressing B-RafV600E.

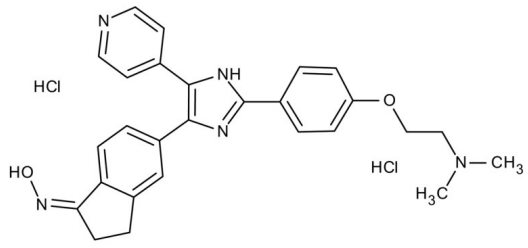
Ordering Information

[Order Online »](#)

ENZ-CHM113-0005	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,N-dimethyl-2-[4-[(4E)-4-(1-nitroso-2,3-dihydroinden-5-ylidene)-5-(1H-pyridin-4-ylidene)-1H-imidazol-2-yl]phenoxy]ethanamine
Appearance	Yellow powder solid.
CAS	405554-55-4
Couple Target	Erk, Raf
Couple Type	Inhibitor
Formula	$C_{27}H_{27}N_5O_2 \cdot HCl$
MW	539
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO.

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-be@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