

Lavendustin A

EGFR inhibitor

Cell permeable tyrosine kinase inhibitor with little effect on cAMP-dependent protein kinase (PKA) or protein kinase C (PKC). Potent and selective inhibitor of the EGF receptor tyrosine kinase ($IC_{50}=11\text{nM}$). Binds to a site on the kinase which is distinct from the ATP and peptide substrate binding sites. Inhibits EGF-induced proliferation of cultured human myometrial smooth muscle cells and suppresses VEGF-induced angiogenesis in rats. Effective at $10\mu\text{M}$ in cultured leech neurons. Also inhibits NMDA-stimulated cGMP production.

Citations: 12

[View Online »](#)

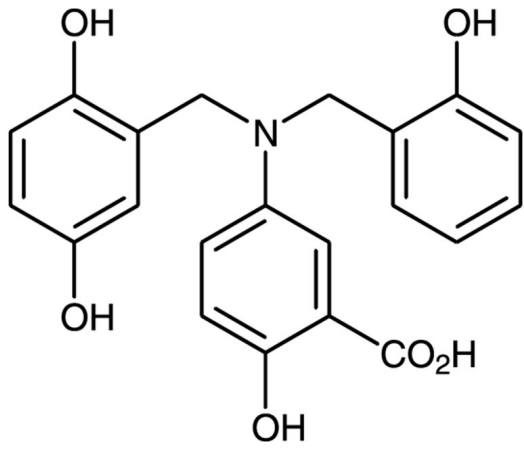
Ordering Information

[Order Online »](#)

BML-EI185-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C.
Handling	Protect from oxygen.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-Amino-[(N-2,5-dihydroxybenzyl)-N'-2-hydroxybenzyl]salicylic acid, RG14355
Appearance	Off-white to tan solid.
CAS	125697-92-9
Couple Target	EGFR
Couple Type	Inhibitor
Formula	$C_{21}H_{19}NO_6$
MW	381.4
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO or ethanol.

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