

SL-327

MEK inhibitor

A selective inhibitor of MEK1 and MEK2 (IC_{50} = 0.18 and 0.22 μ M respectively). Displays neuroprotective effects against ischemic brain injury.

Citations: 7

[View Online »](#)

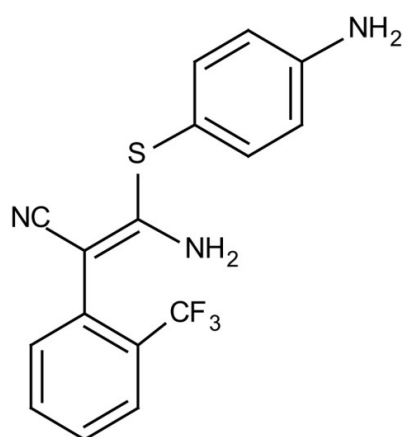
Ordering Information

[Order Online »](#)

BML-EI365-0010	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

Alternative Name [Amino[(4-aminophenyl)thio]methylene]-2-(trifluoromethyl)benzeneacetonitrile

Appearance White to off-white solid.

CAS 305350-87-2

Couple Target MEK

Couple Type Inhibitor

Formula $C_{16}H_{12}F_3N_3S$

MW 335.4

Purity ≥99% (HPLC)

Solubility Soluble in DMSO to 100 mM or in ethanol to 100 mM.

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