

L-744,832

Farnesyltransferase inhibitor

Ras farnesyltransferase inhibitor. Induces tumor regression in transgenic mice with multiple oncogenic mutations by mediating alterations in both cell cycle control and apoptosis. Induces p21 expression and arrests cells at G1. Causes enhanced mitotic sensitivity to taxol.

Citations: 8

[View Online »](#)

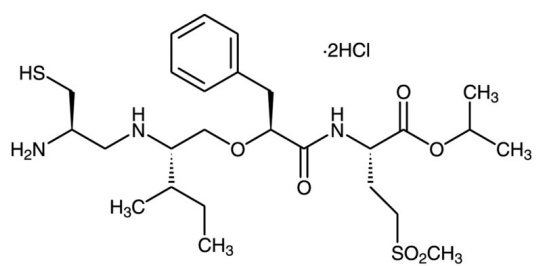
Ordering Information

[Order Online »](#)

BML-G242-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. 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

Appearance	White solid.
CAS	1177806-11-9
Couple Target	Farnesyltransferase
Couple Type	Inhibitor
Formula	$C_{26}H_{47}Cl_2N_3O_6S_2$
MW	632.7
Purity	≥98% (TLC)
Solubility	Soluble in DMSO, water, or acetonitrile.

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-
be@enzolifesciences.com](mailto:info-be@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)