

BI-D1870

S6 kinase inhibitor

Cell permeable, ATP-competitive inhibitor of the four isoforms of ribosomal S6 kinase (p90 RSK) (IC_{50} 's at 100 μ M ATP: RSK1 = 31nM, RSK2 = 24nM, RSK3 = 18nM, RSK4 = 15nM). Does not significantly inhibit 10 other AGC kinases, and over 40 other protein kinases at 100X concentrations. Acts on the N-terminal AGC kinase domain. Completely prevents PMA-induced TSC2 phosphorylation.

Citations: 19

[View Online »](#)

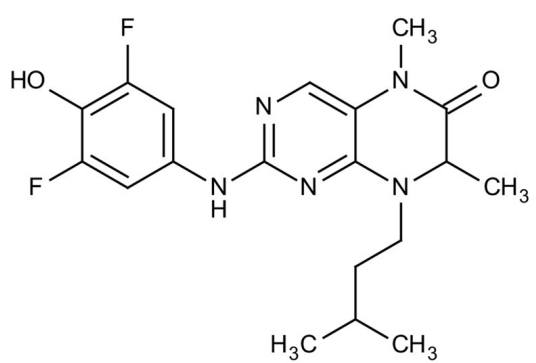
Ordering Information

[Order Online »](#)

BML-EI407-0001	1mg
BML-EI407-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year from the date of receipt when stored as supplied at -20°C
Handling	Solutions can be stored at -20°C for up to 3 months. Protect from moisture and light.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-[(3,5-Difluoro-4-hydroxyphenyl)amino]-7,8-dihydro-5,7-dimethyl-8-(3-methylbutyl)-6(5H)-pteridone
Appearance	Off-white solid.
CAS	501437-28-1
Couple Target	S6 kinase
Couple Type	Inhibitor
Formula	$C_{19}H_{23}F_2N_5O_2$
MW	391.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (>25mg/ml).



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)