

PD0325901

Selective MEK inhibitor

PD0325901 (PD325901) is selective and non ATP-competitive MEK inhibitor with an IC_{50} of 0.33 nM, roughly 500-fold more potent than CI-1040 on the phosphorylation of ERK1 and ERK2.

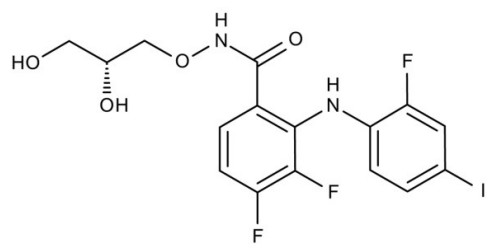
Ordering Information

[Order Online »](#)

ENZ-CHM202-0010	10mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(R)-N-(2,3-Dihydroxypropoxy)-3,4-difluoro-2-((2-fluoro-4-iodophenyl)amino)benzamide
Appearance	White to off-white powder.
CAS	391210-10-9
Couple Target	MEK
Couple Type	Inhibitor
Formula	$C_{16}H_{14}F_3IN_2O_4$
Identity	Determined by NMR.
MW	482.2
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (12mg/ml) or ethanol (6.3mg/ml) with slight warming.

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