

PD0325901

Reprogramming Enhancer -MEK Inhibitor

Catalog # inh-pd32

For research use only

Version # 10C17-MM

PRODUCT INFORMATION

Content:

- 2 mg PD0325901

Storage and stability:

- PD0325901 is provided as a crystalline solid and shipped at room temperature. Store at -20°C. Solid product is stable 1 year at -20°C. Protect from light.
- Upon resuspension in an organic solvent (DMSO/ethanol/DMF), PD0325901 should be aliquoted and stored at -20°C. Avoid repeated freeze-thaw cycles. Resuspended product is stable for 3 months at -20°C when properly stored.
- An aqueous solution can be obtained by diluting the organic solvent into aqueous buffer or isotonic saline. We do not recommend storing the aqueous solution for more than 1 day.

CHEMICAL PROPERTIES

CAS number: 391210-10-9

Synonym: N-[(2R)-2,3-dihydroxypropoxy]-3,4-difluoro-2-[(2-fluoro-4-iodophenyl)amino]-benzamide

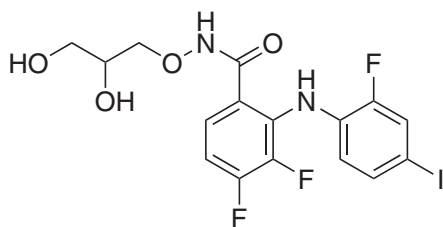
Formula: C₁₆H₁₄F₃IN₂O₄

Molecular weight: 482.19

Solubility: DMSO, ethanol, DMF (20 mg/ml)

Appearance: White to off-white solid

Purity: >98% (HPLC)



DESCRIPTION

PD0325901 is a synthetic organic molecule that selectively binds to and inhibits mitogen-activated protein kinase kinase (MEK). The MEK/ERK signaling pathway plays an important role in the self-renewing state of ES cells¹. Inhibition of this pathway combined with the inhibition of other signaling pathways has been shown to improve the efficiency of iPS cell generation. Combination of PD0325901 with a TGF- β receptor inhibitor was found to improve the reprogramming of OSKM-infected human fibroblasts².

1. Ying QL. *et al.*, 2008. The ground state of embryonic stem cell self-renewal. *Nature*. 453(7194):519-23.

2. Lin T. *et al.*, 2009. A chemical platform for improved induction of human iPSCs. *Nat Methods*. 6(11):805-8.

Working concentration: 0.5 μ M

RELATED PRODUCTS

Product	Catalog Code
5-Aza-cytidine	inh-aza
Bix-01294	inh-bix
SB431542	inh-sb43
Valproic acid	inh-vpa
LENTI-Smart™ OSKM (human)	ltsint-hoskm-5
LENTI-Smart™ OSKM (mouse)	ltsint-moskm-5

TECHNICAL SUPPORT

Toll free (US): 888-457-5873

Outside US: (+1) 858-457-5873

Europe: +33 562-71-69-39

E-mail: info@invivogen.com

Website: www.invivogen.com



3950 Sorrento Valley Blvd. Suite 100
San Diego, CA 92121 - USA