

# SB-202190

## p38 inhibitor

Cell permeable, potent and selective p38 MAP kinase inhibitor ( $IC_{50}=30$  nM). Inhibits p38 $\alpha$  and  $\beta$ , but not  $\gamma$  and  $\delta$  isoforms. Does not inhibit ERK2 or other members of the MAP kinase family or their upstream activators. A useful tool for dissecting the role of p38 in signaling pathways. Induces apoptosis.

Citations: 26

[View Online »](#)

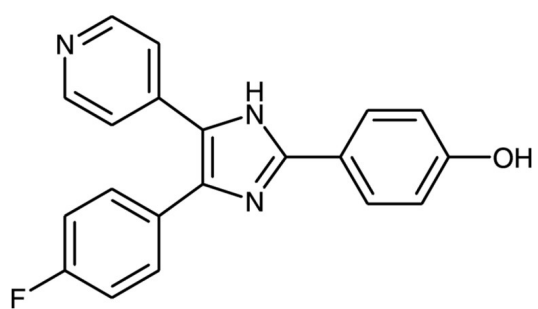
## Ordering Information

[Order Online »](#)

BML-EI294-0001	1mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	FHPI, 4-(4-Fluorophenyl)-2-(4-hydroxyphenyl)-5-(4-pyridyl)-1H-imidazole
Appearance	Pale yellow solid.
CAS	152121-30-7
Couple Target	p38
Couple Type	Inhibitor
Formula	$C_{20}H_{14}FN_3O$
MW	331.3
Purity	≥98% (TLC)
Solubility	Soluble in methanol, DMSO or acetone; slightly soluble in ethyl acetate.

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