

SB202190 . hydrochloride

p38 inhibitor

A water soluble form of the potent p38 (SAPK2a) inhibitor SB202190 (Prod. No. BML-EI294).

Citations: 10

[View Online »](#)

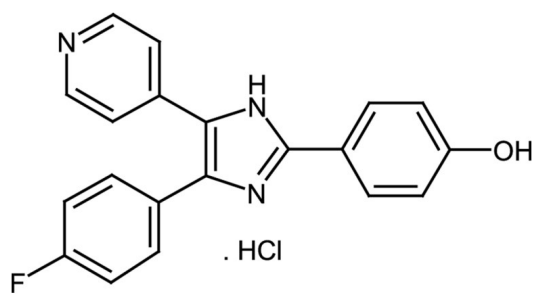
Ordering Information

[Order Online »](#)

ALX-270-268-M001	1mg
ALX-270-268-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	FHPI . HCl, 4-(4-Fluorophenyl)-2-(4-hydroxyphenyl)-5-(4-pyridyl)-1H-imidazole . HCl
Appearance	Yellow solid.
CAS	350228-36-3
Couple Target	p38
Couple Type	Inhibitor
Formula	$C_{20}H_{14}FN_3O \cdot HCl$
Identity	Determined by ¹ H-NMR.
MW	331.3 . 36.5
Purity	≥96% (HPLC)
Solubility	Soluble in water, methanol or DMSO.

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-be@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