

CC-401 . HCl

Potent inhibitor of JNK

CC-401 is a competitive inhibitor of the ATP binding site in the active, phosphorylated, form of JNK. This prevents JNK from phosphorylating its various target molecules, including the amino terminus of c-Jun. It is a potent inhibitor of all three forms of JNK (K_i of 25-50 nM), and has at least 40-fold selectivity for JNK compared with other related kinases, including: p38, ERK, IKK2, PKC, Lck, and ZAP70. CC-401 acts to inhibit JNK signaling by competitive binding to the adenosine triphosphate-binding site in the active, phosphorylated, form of JNK, resulting in inhibition of the phosphorylation of JNK targets, such as the amino-terminal activation domain of the transcription factor, c-Jun. In cell-based assays, 1–5 $\mu\text{mol/l}$ CC-401 provides specific JNK inhibition.

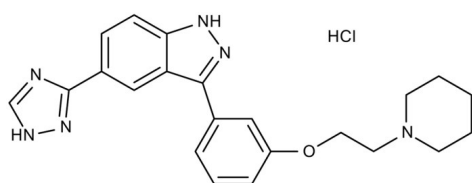
Ordering Information

[Order Online »](#)

ENZ-CHM194-0001	1mg
-----------------	-----

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	3-[3-(2-piperidin-1-ylethoxy)phenyl]-5-(1H-1,2,4-triazol-5-yl)-1H-indazole hydrochloride
Appearance	White powder.
CAS	1438391-30-0
Couple Target	Jnk
Couple Type	Inhibitor
Formula	$C_{22}H_{24}N_6O \cdot HCl$
Identity	Determined by EM-MS, NMR
MW	424.9
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (15mM).



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