

BX795

Inhibitor of PDK1, TBK1, IKK- ϵ , IRF3

BX795 was initially developed as a PDK1 inhibitor (direct competitive inhibitor $IC_{50} = 11\text{nM}$ for PDK1) and has been shown to be a potent and relatively specific inhibitor of TBK1 and IKK- ϵ . It blocks the phosphorylation, nuclear translocation, and transcriptional activity of interferon regulatory factor 3. BX795 also is a potent inhibitor of cell growth of multiple cancer cell lines with IC_{50} values ranging from submicromolar amounts ($0.368\text{ }\mu\text{M}$ to greater than $450\text{ }\mu\text{M}$).

Citations: 4

[View Online »](#)

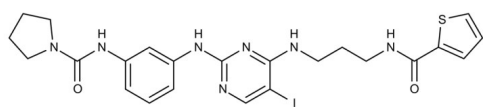
Ordering Information

[Order Online »](#)

ENZ-CHM189-0005	5mg
ENZ-CHM189-0010	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	N-[3-[[5-iodo-4-[3-(thiophene-2-carbonylamino)propylamino]pyrimidin-2-yl]amino]phenyl]pyrrolidine-1-carboxamide
Appearance	White to off-white powder.
CAS	702675-74-9
Couple Target	IKK, IRF, Pyruvate dehydrogenase kinase, TANK-binding kinase
Couple Type	Inhibitor
Formula	$C_{23}H_{26}IN_7O_2S$
Identity	Determined by NMR.
MW	591.5
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (80mg/ml).



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