

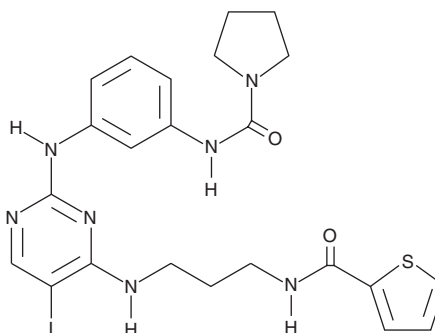
# PRODUCT INFORMATION



## BX-795

Item No. 14932

**CAS Registry No.:** 702675-74-9  
**Formal Name:** N-[3-[[5-iodo-4-[[3-[(2-thienylcarbonyl)amino]propyl]amino]-2-pyrimidinyl]amino]phenyl]-1-pyrrolidinecarboxamide  
**MF:** C<sub>23</sub>H<sub>26</sub>IN<sub>7</sub>O<sub>2</sub>S  
**FW:** 591.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 247 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

BX-795 is supplied as a crystalline solid. A stock solution may be made by dissolving the BX-795 in the solvent of choice, which should be purged with an inert gas. BX-795 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of BX-795 in these solvents is approximately 11.1 and 12.5 mg/ml, respectively.

BX-795 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BX-795 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BX-795 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

BX-795 is a multi-kinase inhibitor that is an ATP-competitive inhibitor of 3-phosphoinositide-dependent protein kinase 1 (PDK1) and TANK-binding kinase 1 (TBK1; IC<sub>50</sub>s = 6 and 11 nM, respectively).<sup>1,2</sup> It also inhibits the CAMK family kinases AMPK-related kinase 5 (NUAK1; IC<sub>50</sub> = 5 nM) and microtubule affinity-regulating kinase 1 (MARK1), MARK2, MARK3, and MARK4 (IC<sub>50</sub>s = 55, 53, 81, and 19 nM, respectively), as well as the MAP3K kinases mixed-lineages kinase 1 (MLK1), MLK2, and MLK3 (IC<sub>50</sub>s = 50, 46, and 42 nM, respectively). BX-795 is selective for these kinases over five additional kinases (IC<sub>50</sub>s = >1,100 nM) but does inhibit the receptor tyrosine kinase VEGFR, as well as IκB kinase ε (IKKε) and Aurora B kinase (IC<sub>50</sub>s = 41 and 31 nM, respectively).<sup>1</sup> It inhibits phosphorylation of Akt at Thr<sup>308</sup> and p70 ribosomal S6 kinase 1 (p70S6K1) at Thr<sup>389</sup> in PC3 prostate cancer cells (IC<sub>50</sub> = 300 nM).<sup>1</sup> BX-795 (1 μM) also inhibits the secretion of IFN-β induced by LPS or poly(I:C) in RAW 264.7 macrophages.<sup>2</sup> It inhibits the proliferation of PC3 prostate and MDA-MB-468 breast cancer cells (IC<sub>50</sub>s = 250 and 720 nM, respectively). BX-795 (5 μM) inhibits the proliferation of PDK1<sup>-/-</sup> murine embryonic stem cells expressing either wild-type PDK1 or PDK1 with an alanine-to-glycine (L159G) mutation.<sup>3</sup>

### References

1. Feldman, R.I., Wu, J.M., Polokoff, M.A., *et al.* Novel small molecule inhibitors of 3-phosphoinositide-dependent kinase-1. *J. Biol. Chem.* **280**(20), 19867-19874 (2005).
2. Clark, K., Plater, L., Pegg, M., *et al.* Use of the pharmacological inhibitor BX795 to study the regulation and physiological roles of TBK1 and IκB kinase ε: A distinct upstream kinase mediates Ser-172 phosphorylation and activation. *J. Biol. Chem.* **284**(21), 14136-14146 (2009).
3. Tamgüney, T., Zhang, C., Fiedler, D., *et al.* Analysis of 3-phosphoinositide-dependent kinase-1 signaling and function in ES cells. *Exp. Cell Res.* **314**(11-12), 2299-2312 (2008).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/04/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM