

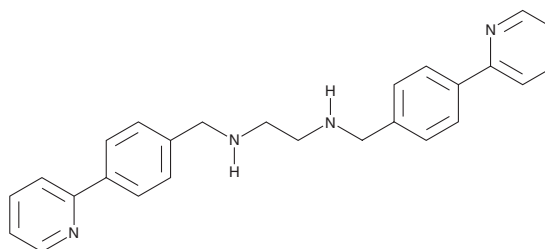
PRODUCT INFORMATION



BC-1215

Item No. 21964

CAS Registry No.: 1507370-20-8
Formal Name: N¹,N²-bis[[4-(2-pyridinyl)phenyl]methyl]-1,2-ethanediamine
MF: C₂₆H₂₆N₄
FW: 394.5
Purity: ≥98%
UV/Vis.: λ_{max}: 250, 280 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

BC-1215 is supplied as a crystalline solid. A stock solution may be made by dissolving the BC-1215 in the solvent of choice. BC-1215 is soluble in the organic solvent dimethyl formamide, which should be purged with an inert gas, at a concentration of approximately 0.5 mg/ml.

Description

BC-1215 is an inhibitor of F-box protein 3 (FBXO3; IC₅₀ = 0.9 µg/mL for Il-1β release).¹ It decreases FBXO3 interaction with F-box/LRR-repeat protein 2 (FBXL2) in a dose-dependent manner and prevents FBXO3-dependent ubiquitination of FBXL2. It also lowers expression of Tnf receptor-associated factors (TRAFs) 1-6 without affecting TRAF1-6 mRNA levels in resting and LPS-stimulated peripheral blood mononuclear cells (PBMCs). BC-1215 (100 µg, i.p., per animal) reduces plasma levels of the inflammatory cytokines Il-6, Tnf, and Il-1β and decreases peritoneal bacteria counts in a mouse model of cecal ligation and puncture-induced sepsis. It also reduces inflammation in mouse models of DSS-induced colitis, paw edema, and ear edema.²

References

1. Chen, B.B., Coon, T.A., Glasser, J.R., *et al.* A combinatorial F box protein directed pathway controls TRAF adaptor stability to regulate inflammation. *Nat. Immunol.* **14**(5), 470-479 (2013).
2. Mallampalli, R.K., Coon, T.A., Glasser, J.R., *et al.* Targeting F box protein Fbxo3 to control cytokine-driven inflammation. *J. Immunol.* **191**(10), 5247-5255 (2013).

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, 11/30/2022

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