

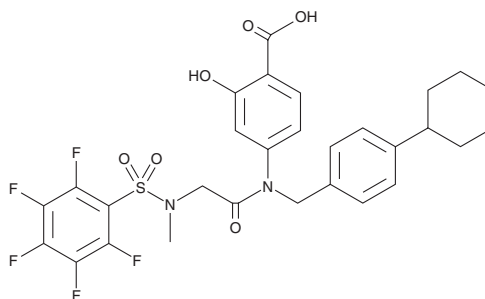
PRODUCT INFORMATION



BP-1-102

Item No. 28368

CAS Registry No.: 1334493-07-0
Formal Name: 4-[[[(4-cyclohexylphenyl)methyl]
[2-[methyl[(2,3,4,5,6-pentafluorophenyl)
sulfonyl]amino]acetyl]amino]-2-hydroxy-
benzoic acid
MF: C₂₉H₂₇F₅N₂O₆S
FW: 626.6
Purity: ≥98%
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

BP-1-102 is supplied as a crystalline solid. A stock solution may be made by dissolving the BP-1-102 in the solvent of choice, which should be purged with an inert gas. BP-1-102 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BP-1-102 in these solvents is approximately 15 and 10 mg/ml, respectively.

Description

BP-1-102 is an inhibitor of STAT3.¹ It binds to STAT3 ($K_d = 504$ nM) and inhibits STAT3 binding to an IL-6R/gp130 peptide in a fluorescence polarization assay ($IC_{50} = 4.1$ μM). BP-1-102 (10-30 μM) decreases viability and proliferation of STAT3-dependent NIH3T3/v-Src, MDA-MB-231, PANC-1, DU145 and A549 cancer cells, but not STAT3-independent NIH3T3, NIH3T3/vRas, TE-71, and A2780S cancer cells. It reduces tumor growth and inhibits expression of the STAT3-dependent genes encoding c-Myc, survivin, Bcl-xL, cyclin D1, and VEGF in MDA-MB-231 and A549 mouse xenograft models when administered at doses of 1 and 3 mg/kg.

Reference

1. Zhang, X., Yue, P., Page, B.D., *et al.* Orally bioavailable small-molecule inhibitor of transcription factor Stat3 regresses human breast and lung cancer xenografts. *Proc. Natl. Acad. Sci. USA.* **109(24)**, 9623-9628 (2012).

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, 11/07/2022

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