

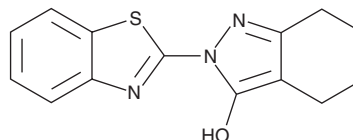
PRODUCT INFORMATION



BD 750

Item No. 37574

CAS Registry No.: 892686-59-8
Formal Name: 2-(2-benzothiazolyl)-4,5,6,7-tetrahydro-2H-indazol-3-ol
MF: C₁₄H₁₃N₃OS
FW: 271.3
Purity: ≥98%
Supplied as: A 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

BD 750 is supplied as a solid. A stock solution may be made by dissolving the BD 750 in the solvent of choice, which should be purged with an inert gas. BD 750 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BD 750 in DMSO and DMF is approximately 30 and 10 mg/ml, respectively. BD 750 is slightly soluble in ethanol.

Description

BD 750 is an immunosuppressant.¹ It inhibits the proliferation of primary human and mouse T cells (IC_{50} s = 1.1 and 1.5 μ M, respectively) without inducing cytotoxicity when used at concentrations ranging from 3.13 to 50 μ M. BD 750 induces cell cycle arrest at the G₀/G₁ phases in CD3 and CD28 antibody-activated primary human T cells when used at concentrations of 0.625, 2.5, and 10 μ M and decreases the levels of cyclin D3 and cyclin-dependent kinase 6 (Cdk6) in CD3 and CD28 antibody-activated primary mouse T cells at 2.5 and 10 μ M. It reduces the proliferation of IL-2-stimulated CTLL mouse T cells (IC_{50} = 1.2 μ M) and decreases phosphorylated Stat5 levels in the same cells when used at concentrations of 5 and 20 μ M. *In vivo*, BD 750 inhibits increases in ear thickness and ear patch weight in a mouse model of delayed-type hypersensitivity induced by 2,4-dinitrofluorobenzene.

Reference

1. Liu, Y., Yang, T., Li, H., *et al.* BD750, a benzothiazole derivative, inhibits T cell proliferation by affecting the JAK3/STAT5 signalling pathway. *Br. J. Pharmacol.* **168**(3), 632-643 (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, 10/30/2023

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