

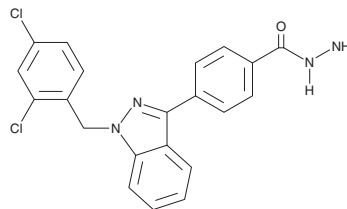
# PRODUCT INFORMATION



## BHD

Item No. 38962

**CAS Registry No.:** 2749517-98-2  
**Formal Name:** 4-[1-[(2,4-dichlorophenyl)methyl]-1H-indazol-3-yl]-benzoic acid hydrazide  
**MF:** C<sub>21</sub>H<sub>16</sub>Cl<sub>2</sub>N<sub>4</sub>O  
**FW:** 411.3  
**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

BHD is supplied as a solid. A stock solution may be made by dissolving the BHD in the solvent of choice, which should be purged with an inert gas. BHD is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BHD in these solvents is approximately 10 and 20 mg/ml, respectively.

### Description

BHD is a reversible male contraceptive agent.<sup>1</sup> *In vivo*, BHD (100 and 500 mg/kg) induces spermatogenic cell apoptosis and seminiferous tubule abnormalities, disrupts the blood-testis barrier, and reduces the fertility index in male mice for a duration of two to six weeks post-administration.

### Reference

1. Gong, S., Zhu, S., Zhou, P., et al. Discovering a reversible male contraceptive agent derived from Ionidamine. *ACS Omega* **8**(20), 18245-18254 (2023).

#### 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, 09/12/2025

#### 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](http://WWW.CAYMANCHEM.COM)