

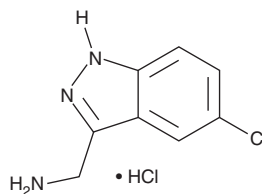
PRODUCT INFORMATION



BI SF3 (hydrochloride)

Item No. 41113

CAS Registry No.: 1801257-24-8
Formal Name: 5-chloro-1H-indazole-3-methanamine, monohydrochloride
MF: $C_8H_8ClN_3 \cdot HCl$
FW: 218.1
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

BI SF3 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the BI SF3 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. BI SF3 (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of BI SF3 (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. BI SF3 (hydrochloride) is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

BI SF3 is a synthetic intermediate.¹ It has been used in the synthesis of quinolinylloxazoles with kallikrein inhibitory activity.

Reference

1. McDonald, A. and Qian, S. Therapeutic inhibitory compounds. *Lifesci Pharmaceuticals, Inc.* WO2018/011628A1 (2018).

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, 09/24/2024

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