

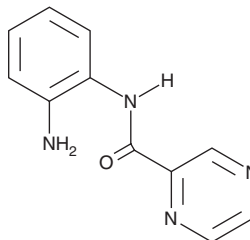
PRODUCT INFORMATION



BG45

Item No. 16918

CAS Registry No.: 926259-99-6
Formal Name: N-(2-aminophenyl)-2-pyrazinecarboxamide
MF: C₁₁H₁₀N₄O
FW: 214.2
Purity: ≥95%
UV/Vis.: λ_{max}: 267 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

BG45 is supplied as a crystalline solid. A stock solution may be made by dissolving the BG45 in the solvent of choice, which should be purged with an inert gas. BG45 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BG45 in these solvents is approximately 20, 30, and 5 mg/ml, respectively.

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 BG45 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of BG45 in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

BG45 is an HDAC class I inhibitor with selectivity for HDAC3 (IC₅₀ = 289 nM). It inhibits HDAC1, HDAC2, and HDAC6 with greatly reduced potency (IC₅₀s = 2, 2.2, and >20 μM, respectively).¹ At concentrations up to 50 mg/kg, BG45 alone or in combination with bortezomib has been shown to dose-dependently inhibit tumor growth in a mouse model of multiple myeloma.¹

Reference

1. Minami, J., Suzuki, R., Mazitschek, R., *et al.* Histone deacetylase 3 (HDAC3) as a novel therapeutic target in multiple myeloma. *Leukemia* **28**(3), 680-689 (2014).

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, 11/23/2022

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