

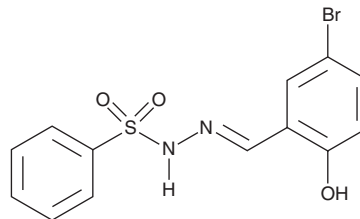
PRODUCT INFORMATION



Shz-1

Item No. 15211

CAS Registry No.: 326886-05-9
Formal Name: benzenesulfonic acid, 2-[(5-bromo-2-hydroxyphenyl)methylene]hydrazide
Synonym: Sulfonylhydrazone-1
MF: C₁₃H₁₁BrN₂O₃S
FW: 355.2
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 275, 329 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

Shz-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the Shz-1 in the solvent of choice, which should be purged with an inert gas. Shz-1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Shz-1 in these solvents is approximately 30 mg/ml.

Description

Shz-1 is a Shz that activates cardiac differentiation, inducing the expression of early cardiac genes in mouse pluripotent and embryonic stem cells *in vitro* at low micromolar concentrations.¹ In addition, Shz-1 drives differentiation of human peripheral blood mononuclear cells, mobilized by granule colony-stimulating factor, toward a cardiac phenotype.¹ These Shz-1 treated cells enhance myocardial regenerative repair in cryo-injured rat heart.¹

Reference

1. Sadek, H., Hannack, B., Choe, E., *et al.* Cardiogenic small molecules that enhance myocardial repair by stem cells. *Proc. Natl. Acad. Sci. USA* **105(16)**, 6063-6068 (2008).

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, 10/19/2022

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