

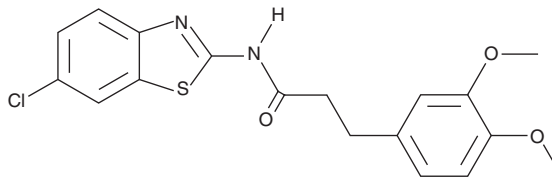
PRODUCT INFORMATION



KY 02111

Item No. 14315

CAS Registry No.: 1118807-13-8
Formal Name: N-(6-chloro-2-benzothiazolyl)-3,4-dimethoxy-benzenepropanamide
MF: C₁₈H₁₇ClN₂O₃S
FW: 376.9
Purity: ≥95%
UV/Vis.: λ_{max}: 204, 223, 281, 304 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

KY 02111 is supplied as a crystalline solid. A stock solution may be made by dissolving the KY 02111 in the solvent of choice, which should be purged with an inert gas. KY 02111 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

KY 02111 promotes the differentiation of human pluripotent stem cells to cardiomyocytes by inhibiting Wnt signaling.¹ Treatment with 10 μM of KY 02111 has been shown to increase the ratio of beating cardiac colonies by 70–94% in cell aggregates of the embryonic stem cell lines KhES-1 and KhES-3 as well as the induced pluripotent stem cell lines 253G1, IMR90-1, IMR90-4, and RCHIPC0003.¹

Reference

1. Minami, I., Yamada, K., Otsuji, T.G., *et al.* A small molecule that promotes cardiac differentiation of human pluripotent stem cells under defined, cytokine- and xeno-free conditions. *Cell Rep.* **2**(5), 1448-1460 (2012).

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, 06/29/2020

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