

PRODUCT INFORMATION



KC159

Item No. 43640

Formal Name: 4-(5-((2,6-dichlorobenzyl)thio)-1,3,4-thiadiazol-2-yl)benzoic acid

MF: C₁₆H₁₀Cl₂N₂O₂S₂

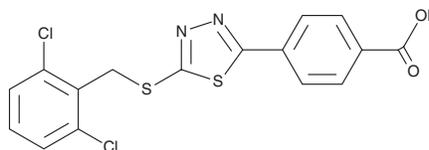
FW: 397.3

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

KC159 is supplied as a solid. A stock solution may be made by dissolving the KC159 in the solvent of choice, which should be purged with an inert gas. KC159 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

KC159 is an agonist of Piezo1.¹ It induces calcium influx in HEK cells overexpressing mouse Piezo1 (EC₅₀ = 280 nM) and induces inward currents in a whole-cell patch-clamp assay using HEK cells overexpressing mouse Piezo1 when used at a concentration of 10 μM. KC159 induces smooth muscle relaxation in isolated and perfused mouse portal veins (EC₅₀ = 1.14 μM).

Reference

1. Parsonage, G., Cuthbertson, K., Endesh, N., *et al.* Improved PIEZO1 agonism through 4-benzoic acid modification of Yoda1. *Br. J. Pharmacol.* **180(16)**, 2039-2063 (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, 06/02/2025

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