

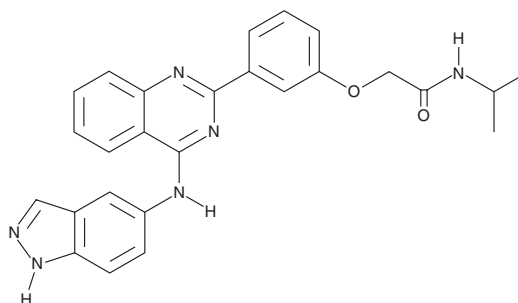
# PRODUCT INFORMATION



## KD 025

Item No. 17055

**CAS Registry No.:** 911417-87-3  
**Formal Name:** 2-[3-[4-(1H-indazol-5-ylamino)-2-quinazoliny]phenoxy]-N-(1-methylethyl)-acetamide  
**Synonyms:** Belumosudil, SLx-2119  
**MF:** C<sub>26</sub>H<sub>24</sub>N<sub>6</sub>O<sub>2</sub>  
**FW:** 452.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 248, 304, 338 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

KD 025 is supplied as a crystalline solid. A stock solution may be made by dissolving the KD 025 in the solvent of choice, which should be purged with an inert gas. KD 025 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of KD 025 in these solvents is approximately 25 and 30 mg/ml, respectively.

KD 025 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

## Description

KD 025 is an inhibitor of Rho-associated kinase 2 (ROCK2; IC<sub>50</sub> = 0.105 μM).<sup>1</sup> It is selective for ROCK2 over ROCK1 (IC<sub>50</sub> = 24 μM). KD 025 (10 μM) decreases the expression of connective tissue growth factor (CTGF) and induces remodeling of the actin cytoskeleton in isolated human ileal fibrotic smooth muscle cells. It inhibits heat-killed *C. albicans*- or *S. epidermidis*-induced production of IL-17 in isolated human peripheral blood mononuclear cells (PBMCs) when used at concentrations ranging from 1.25 to 10 μM.<sup>2</sup> KD 025 (100 and 200 mg/kg) reduces infarct volume in a mouse model of cerebral ischemia induced by transient middle cerebral artery occlusion (MCAO).<sup>4</sup> It decreases disease severity in a mouse model of sclerodermatous chronic graft versus host disease (GVHD) when administered at a dose of 150 mg/kg.<sup>3</sup> Formulations containing KD 025 have been used in the treatment of chronic GVHD.

## References

1. Boerma, M., Fu, Q., Wang, J., *et al.* Comparative gene expression profiling in three primary human cell lines after treatment with a novel inhibitor of Rho kinase or atorvastatin. *Blood Coagul. Fibrinolysis* **19**(7), 709-718 (2008).
2. Tengesdal, I.W., Kitzenberg, D., Li, S., *et al.* The selective ROCK2 inhibitor KD025 reduces IL-17 secretion in human peripheral blood mononuclear cells independent of IL-1 and IL-6. *Eur. J. Immunol.* **48**(10), 1679-1686 (2018).
3. Flynn, R., Paz, K., Du, J., *et al.* Targeted Rho-associated kinase 2 inhibition suppresses murine and human chronic GVHD through a Stat3-dependent mechanism. *Blood* **127**(17), 2144-2154 (2016).
4. Wang, H., Zhang, D., Ge, M., *et al.* Formononetin inhibits enterovirus 71 replication by regulating COX-2/PGE<sub>2</sub> expression. *Virology* **12**, 35 (2015).

### 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, 10/26/2022

## 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