

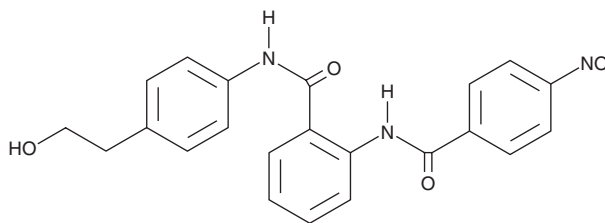
PRODUCT INFORMATION



KS-176

Item No. 22201

CAS Registry No.: 1253452-78-6
Formal Name: N-[4-(2-hydroxyethyl)phenyl]-
2-[(4-nitrobenzoyl)amino]-
benzamide
MF: C₂₂H₁₉N₃O₅
FW: 405.4
Purity: ≥98%
UV/Vis.: λ_{max}: 260 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

KS-176 is supplied as a crystalline solid. A stock solution may be made by dissolving the KS-176 in the solvent of choice, which should be purged with an inert gas. KS-176 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of KS-176 in these solvents is approximately 20 mg/ml.

KS-176 is sparingly soluble in aqueous buffers. KS-176 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

KS-176 is a selective inhibitor of breast cancer resistance protein (BCRP; IC₅₀ = 1.39 μM) with no activity against P-glycoprotein (P-gp) or multidrug resistance-associated protein 1 (MRP1).¹ It acts as a non-competitive inhibitor whose potency is increased by imatinib binding. KS-176 does not reduce cell viability at concentrations up to 10 μM.

Reference

1. Pick, A., Klinkhammer, W., and Wiese, M. Specific inhibitors of the breast cancer resistance protein (BCRP). *ChemMedChem* **5**(9), 1498-1505 (2010).

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