

KS-176

BCRP inhibitor

A potent specific inhibitor of Breast Cancer Resistance Protein (BCRP) / ATP-binding cassette sub-family G member 2 (ABCG2).

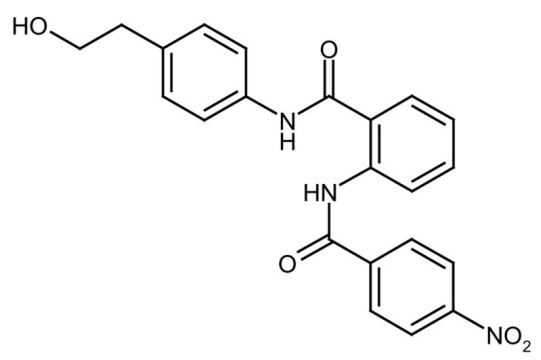
IC_{50} = 1.39 μ M

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -20°C. Stock solutions are stable for up to 2 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(4-(2-hydroxyethyl)phenyl)-2-(4-nitrobenzamido)benzamide
Appearance	Pale yellow solid.
Couple Target	BCRP
Couple Type	Inhibitor
Formula	$C_{22}H_{19}N_3O_5$
MW	405.4
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol, methanol or DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com