

Ko143

BCRP inhibitor

Analog of the fungal toxin fumitremorgin C (Prod. No. ALX-350-127). Potent and selective inhibitor of the breast cancer resistance protein multidrug transporter (BCRP, $EC_{50}=26\text{nM}$). Highly effective for increasing the intracellular drug accumulation and reversing BCRP-mediated multidrug resistance yet has little activity against P-glycoprotein, multidrug resistance-associated protein (MRP1) and other known drug transporters. In mouse models Ko143 was able to restore the intracellular accumulation and abolish the directionality in net flux of dasatinib. It is a highly useful tool for the identification of drug transport processes in complex experimental systems.

Citations: 10

[View Online »](#)

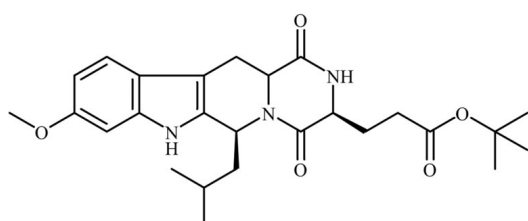
Ordering Information

[Order Online »](#)

BML-EI396-0005	5mg
BML-EI396-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(3S,6S,12aS)-1,2,3,4,6,7,12,12a-Octahydro-9-methoxy-6-(2-methylpropyl)-1,4-dioxopyrazino-[1',2':1,6]pyrido[3,4-b]indole-3-propanoic acid 1,1-dimethylethyl ester
Appearance	White solid.
CAS	461054-93-3
Couple Target	BCRP
Couple Type	Inhibitor
Formula	$C_{26}H_{35}N_3O_5$
MW	469.6
Purity	≥96% (HPLC)
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/ml).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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