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, EC₉₀=26nM). 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 »

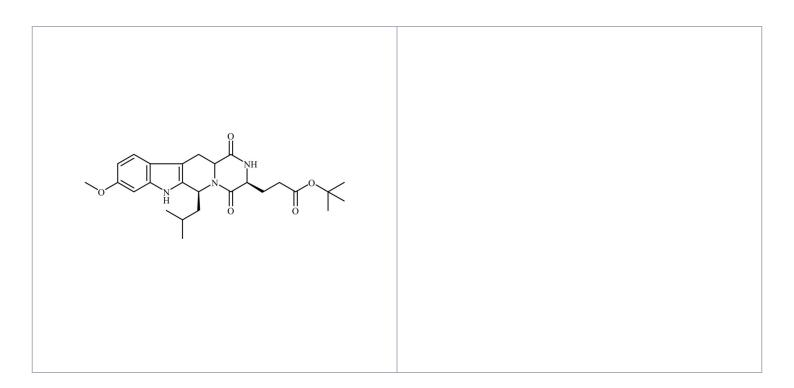
Ordering Information

Order Online »

BML-El396-0005	5mg
BML-El396-0025	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).

Last modified: May 29, 2024

