

Quizartinib

Inhibitor of FLT3 and ABC proteins

Quizartinib (AC220) is a second-generation FLT3 inhibitor that has shown promising activity in Phase II clinical trials. It inhibits mutant and wild-type FLT3 in vivo at 0.1 and 0.5 μM , respectively, and has shown favorable activity and tolerability in phase I and II trials in acute myeloid leukemia, with QT prolongation as the dose-limiting toxicity. Recently, AC220 has also been shown to be an effective inhibitor of ATP-binding cassette (ABC) proteins ABCB1 (P-glycoprotein) and ABCG2 (breast cancer resistance protein). Quizartinib inhibited transport of fluorescent ABCG2 and ABCB1 substrates in ABCG2- and ABCB1-overexpressing cells in a concentration-dependent manner, from 0.1 to 5 μM and from 0.5 to 10 μM , respectively, and inhibited [125I]-IAAP photolabeling of ABCG2 and ABCB1 with IC_{50} values of 0.07 and 3.3 μM , respectively.

Citations: 1

[View Online »](#)

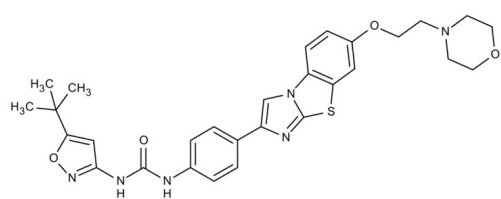
Ordering Information

[Order Online »](#)

ENZ-CHM154-0001	1mg
ENZ-CHM154-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	AC-220, 1-(5-tert-Butyl-1,2-oxazol-3-yl)-3-[4-[6-(2-morpholin-4-ylethoxy)imidazo[2,1-b][1,3]benzothiazol-2-yl]phenyl]urea
Appearance	White powder.
CAS	950769-58-1
Couple Target	BCRP, Flt, P-glycoprotein
Couple Type	Inhibitor
Formula	$C_{29}H_{32}N_6O_4S$
Identity	Determined by EM-MS, NMR
MW	560.7
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (33mg/ml) or ethanol (0.5mg/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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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

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