

E-4031 . dihydrochloride

K⁺ channel blocker

Potent and selective blocker of the inactivating inward rectifying potassium current IKR (HERG channel). Class III antiarrhythmic agent with differential atrial vs. ventricular activity.

Citations: 27

[View Online »](#)

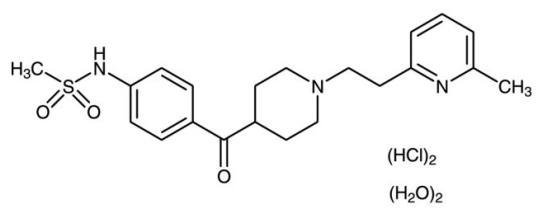
Ordering Information

[Order Online »](#)

BML-KC158-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4'-[[1-[2-(6-methyl-2-pyridyl)ethyl]-4-piperidinyl]carbonyl]-methanesulfonanilide
Appearance	White solid.
CAS	113559-13-0
Couple Target	HERG, Potassium channel
Couple Type	Blocker
Formula	$C_{21}H_{27}N_3O_3S \cdot 2HCl \cdot 2H_2O$
MW	401.5 . 73.0 . 36.0
Purity	≥98% (HPLC)
Solubility	Soluble in any aqueous buffer. Centrifuge all product preparations before use (10000 x g for 5 min).

Last modified: May 29, 2024



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-
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)