

Clofilium . tosylate

K⁺ channel blocker

K⁺ channel blocker. Antiarrhythmic. Cardiac depressant.

Citations: 3

[View Online »](#)

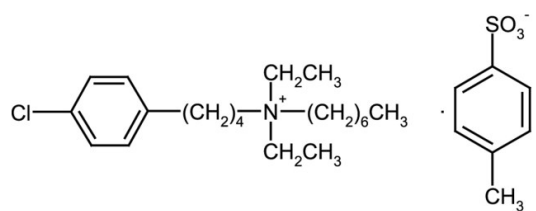
Ordering Information

[Order Online »](#)

ALX-550-242-M100	100mg
ALX-550-242-M250	250mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 4-Chloro-N,N-diethyl-N-heptylbenzenebutanaminium . tosylate

Appearance White solid.

CAS 92953-10-1

Couple Target Potassium channel

Couple Type Blocker

Formula $C_{21}H_{37}ClN \cdot C_7H_7O_3S$

Identity Determined by 1H -NMR.

MW 339.0 . 171.2

Purity $\geq 95\%$ (HPLC)

Solubility Soluble in water, methanol or DMSO.

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)