

Bumetanide

Na⁺ channel blocker

Bumetanide is a specific inhibitor of Na-K-2Cl cotransport. It binds specifically, saturably and with high affinity to membranes from a large variety of cells. A clinically useful diuretic.

Citations: 4

[View Online »](#)

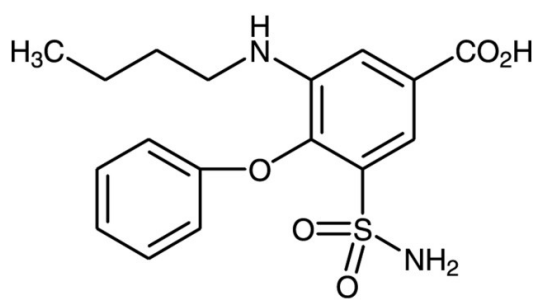
Ordering Information

[Order Online »](#)

BML-EI168-0005	5g
----------------	----

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

Appearance White solid.

CAS 28395-03-1

Couple Target Sodium channel

Couple Type Blocker

Formula $C_{17}H_{20}N_2O_5S$

MW 364.4

Purity $\geq 98\%$

Solubility Soluble in DMSO (25mg/ml) or 100% ethanol (10mg/ml).

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-
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)