

SCREEN-WELL® Ion Channel ligand library

The SCREEN-WELL® Ion Channel Ligand Library contains 70 ion channel blockers and openers for use in characterizing and identifying ion channels in individual cells or tissue.

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|----------------------|------------------------|
| BML-2805-0100 | 100 µl/well - 1Library |
|----------------------|------------------------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|--------------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Stable for at least one year from the date of receipt when stored at -80°C. |
| Long Term Storage | -80°C |
| Shipping | Dry Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|----------------------|---------------|
| Application | HTS |
| Concentration | 10mM in DMSO. |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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