



SCREEN-WELL®

Natural Product library

Natural products are an unsurpassed source of chemical diversity and an ideal starting point for any screening program for pharmacologically active small molecules. Historically, natural products have been the most successful source of new drugs. The SCREEN-WELL® Natural Product library offers over 500 isolated natural compounds of known structure and includes many compounds from ethnopharmacology. Compounds are provided in solution in DMSO. Individual compounds or subsets can be resupplied in gram quantities.

Citations: 28

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|---------------|------------------------|
| BML-2865-0500 | 500 µl/well - 1Library |
| BML-2865-0100 | 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 | 2mg/ml in DMSO |
| Contents | 502 compounds including compounds from the following classes of products: Terpenoids, Peptolides, Flavones, Coumarins, Alkaloids, Macrolides, Isoflavones, Synthetic derivates and many more. |
| Quantity | 500µl/well |

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