



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



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)