

SCREEN-WELL[®] FDA approved drug library V2

Accelerates drug optimization with a unique collection of compounds of known safety

Drug repositioning or repurposing can be an important part of any drug discovery program and has led to several blockbuster drugs (e.g. Viagra and Rogaine). High-content screens, new biomarkers and noninvasive imaging techniques have created new opportunities for pursuing novel indications for approved compounds. Being FDA approved and in use in the clinics, all of the compounds in SCREEN-WELL[®] FDA v. 2.0 Approved Drug Library have known and well-characterized bioactivity, safety and bioavailability - properties which could dramatically accelerate drug development and optimization. The library also avoids irrelevant compounds found in similar libraries, such as herbicides, insecticides, sunscreen agents, and cytotoxic agents. This set provides one of the richest sources of bioactives and offers the greatest degree of drug-likeness available. Hits from this set will provide a significant head start in any drug optimization program. Many of the compounds included can also be provided as standalone compounds for further testing.

Citations: 80

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2843-0100

100 µl/well - 1Library

Manuals, SDS & CofA

[View Online »](#)

- Compounds can be used for diverse drug discovery in the fields of cardiology, neuropsychiatry, immunology, oncology and more!
- Contains more than 100 additional compounds compared to our previous FDA approved library versions
- 100% known bioactivity and greatest degree of drug-likeness available
- Recently approved FDA compounds are also included
- We provide detailed information in our comprehensive documentation for each compound

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 | 10 mM* in DMSO***Compounds are predominantly at concentration of 10mM unless otherwise stated.**Compounds are predominantly provided in DMSO. Selected compounds are dissolved in nuclease-free water. |
| Contents | Over 770 compounds, selected to maximize chemical and pharmacological diversity |
| Quantity | 100µl per 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)