

# SCREEN-WELL®

## Orphan ligand library

The SCREEN-WELL® Orphan ligand library contains 84 compounds with defined, putative, potential or speculative biological activity but whose protein binding partners have not been identified. Ligands are supplied dissolved in biocompatible solvents at 10mM and aliquoted into a deep-well plate at 500µl per well. The library is a rich source of ligands for receptor de-orphaning. Includes the following classes of ligands: Trace Amines, Neurotransmitter Metabolites, Endogenous β-carbolines, Urinary Metabolites, Nicotine Congeners, D-Amino Acids. Receptors for endogenous trace amines have been deorphanized. Endogenous β-carbolines have been shown to display varied biological effects such as hallucinations, hypotension and others. Nicotine congeners such as normicotine are biologically active and may act at novel nicotinic receptor subtypes.

Citations: 9

[View Online »](#)

### Ordering Information

[Order Online »](#)

BML-2825-0500

500 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
Quantity	84 compounds, 500µl per compound

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)