

# SCREEN-WELL<sup>®</sup> ICCB

## Known Bioactives library

The ICCB Known Bioactives Library is an important product for use in chemical genetics and drug discovery. The library is a collection of 472 diverse biologically active compounds with defined biological activity and was developed in collaboration with the Harvard Institute of Chemistry and Cell Biology. It includes the following classes of compounds: GPCR ligands, Second messenger modulators, Nuclear receptor ligands, Actin & tubulin modulators, Kinase inhibitors, Protease inhibitors, Ion channel blockers, Gene regulation agents, Lipid biosynthesis inhibitors, many other classes. Each compound is supplied dissolved in DMSO at 100µL and is ready for assaying. The library can be used for assay development and validation, mechanism profiling, lead screening or as a reference library.

**Citations:** 32

[View Online »](#)

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Application</b>	HTS
<b>Concentration</b>	DMSO solutions (concentration vary by compound and are provided in the documentation)
<b>Contents</b>	472 compounds. Includes compounds that affect most cellular processes and drug target classes, including: GPCR ligands, Second messenger modulators, Nuclear receptor ligands, Actin and tubulin modulators, Kinase inhibitors, Protease inhibitors, Ion channel blockers, Gene regulation agents, Lipid biosynthesis inhibitors, Phosphodiesterase inhibitors, and many other classes.
<b>Quantity</b>	100µl per well

**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)