SCREEN-WELL® 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 »

Ordering Information

Order Online »

BML-2840-0100

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 DMSO solutions (concentration vary by compound and are

provided in the documentation)

Contents 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.

Quantity 100μl per well

Last modified: May 29, 2024

