

# PRODUCT DATA SHEET



## ICCB Known Bioactives Library

**CATALOG NO.:** 2840

**Version:** 2.0

**Lot #:** TEMP

**DESCRIPTION:** BIOMOL has created a novel library of biologically active small organic molecules in cooperation with the Harvard Institute of Chemistry and Cell Biology (ICCB; now the Broad Institute Chemical Biology Program (BCBP) and ICCB-Longwood). Known active compounds are indispensable tools for target validation, biological mechanism profiling and assay validation. The ICCB Known Bioactives Library serves as a widely used standard set of actives for high-throughput screening.

**CONTENTS:** 480 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

*The complete list of compounds is available within the documentation included with the library or by request.*

**FORMAT:** DMSO solutions (concentrations vary by compound and are provided in the documentation). 6 x 96-well plate. 100 µl per well.

**SUPPORTING INFORMATION:** Documentation containing a complete list of compounds and supporting information is supplied with the library or by request at [compoundlibraries@enzolifesciences.com](mailto:compoundlibraries@enzolifesciences.com). Documentation includes physical information, plate positions, and a structure database (SD) file for integration into your compound database.

**STORAGE:** -80°C. Thaw the plates in a 37°C bath.

### REFERENCES & CITATIONS:

1) A. Shelat and K.G. Guy *Nat. Chem. Biol.* 2007 **3** 442

**FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.**