

## Screen-Well™ FDA approved drug library (10 mM)

**CATALOG NO.:** BML-2842

**Version:** 1.3

**Lot #:** TEMP

**DESCRIPTION:** 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 BIOMOL's FDA Approved Drug Library have known and well-characterized bioactivities and have undergone extensive safety & bioavailability testing, which could dramatically accelerate your drug discovery program. 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.

**CONTENTS:** 640 compounds, selected to maximize chemical and pharmacological diversity

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

**FORMAT:** 10 mM DMSO solutions. 8 x 96-well plates. 100µl or 500µl per well.

**SUPPORTING INFORMATION:** Documentation containing a complete list of compounds and supporting information is supplied with the library or by request at [techserv-usa@enzolifesciences.com](mailto:techserv-usa@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.

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