# SCREEN-WELL<sup>®</sup> FDA approved drug library V2

## Accelerates drug optimization with a unique collection of compounds of known safety

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 SCREEN-WELL® FDA v. 2.0 Approved Drug Library have known and well-characterized bioactivity, safety and bioavailability - properties which could dramatically accelerate drug development and optimization. The library also avoids irrelevant compounds found in similar libraries, such as herbicides, insecticides, sunscreen agents, and cytotoxic agents. 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. Many of the compounds included can also be provided as standalone compounds for further testing.

Citations: 80

View Online »

**Ordering Information** 

**Order Online** »

BML-2843-0100

100 µl/well - 1Library

Manuals, SDS & CofA

View Online »

- Compounds can be used for diverse drug discovery in the fields of cardiology, neuropsychiatry, immunology, oncology and more!
- Contains more than 100 additional compounds compared to our previous FDA approved library versions
- 100% known bioactivity and greatest degree of drug-likeness available
- Recently approved FDA compounds are also included
- We provide detailed information in our comprehensive documentation for each compound

### **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 10 mM\* in DMSO\*\*\*Compounds are predominantly at

concentration of 10mM unless otherwise

stated.\*\*Compounds are predominantly provided in DMSO. Selected compounds are dissolved in nuclease-

free water.

Contents Over 770 compounds, selected to maximize chemical and

pharmacological diversity

Quantity 100µl per well

Last modified: May 29, 2024

