# SCREEN-WELL® Cardiotoxicity library

Essential standards for predictive toxicology screening

The SCREEN-WELL<sup>®</sup> Cardiotoxicity Library is a focused collection of 130 compounds with defined and diverse cardiotoxicity, including ion channel blockage, mitochondrial toxicity, arrhythmia, fibrosis, and many more. A variety of structurally and mechanistically different compound classes are included, as well as nontoxic controls. The compounds are dissolved in DMSO at 10mM and aliquoted into deep-well plates. The library is an essential tool for predictive toxicology screening and assay development.

Citations: 6

View Online »

## **Ordering Information**

**Order Online** »

BML-2850-0100	100 μl/well - 1Library
BML-2850-0500	500 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.

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -80°C

**Shipping** Dry Ice

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Application HTS

Concentration DMSO solutions (10mM).

Contents 130 compounds with defined and diverse cardiotoxicity. A

variety of structurally and mechanistically different compound classes, as well as nontoxic controls, are

included.

**Quantity** 100 μl/well or 500 μl/well

Technical Info / Product Notes Application Note (Cardiotoxicity Application):

Multi Parameter in vitro Assessment of Compound Effects on Cardiomyocyte Physiology Using Induced Pluripotent

Stem Cells (iPSC)

Last modified: May 29, 2024

