

PRODUCT DATA SHEET



Screen-Well™ Natural Product Library

CATALOG NO.: 2865

Version: 5.5

Lot #: TEMP

DESCRIPTION: Natural products are an unsurpassed source of chemical diversity and are an ideal starting point for any screening program for pharmacologically active small molecules. Historically, natural products have been the most successful source of new drugs. The Screen-Well™ Natural Product Library includes compounds from microbial, plant and marine sources including many compounds identified from ethnopharmacology (traditional medicine). All of the compounds are of low molecular weight and known structure. The compounds are highly purified. No extracts or mixtures are included in this library.

CONTENTS: 502 compounds including compounds from the following classes of natural products:

Terpenoids	Peptolides	Flavones
Coumarins	Alkaloids	Macrolides
Isoflavones	Synthetic derivatives	
...and many more		

The complete list of compounds is available within the documentation included with the library.

FORMAT: 2 mg/ml DMSO solutions. 6 x 96-well plates. 500µl 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. 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.