

PRODUCT DATA SHEET



Screen-Well™ Rare Natural Product Library

CATALOG NO.: 2866

Version: 1.1

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™ Rare Natural Product Library includes compounds from microbial, plant and marine sources whose supply is limited. 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: 84 compounds

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

Please note: because the compounds in this library are of limited supply and the raw materials can be difficult to obtain, the contents of this library are subject to change frequently, and resupply of individual compounds cannot be guaranteed.

FORMAT: 1 mg/ml DMSO solutions. 1 x 96-well plates. 100µl 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.