

PRODUCT DATA SHEET



Screen-Well™ Orphan Ligand Library

CATALOG NO.: 2825

Version: 3.1

Lot #: SAMPLE

DESCRIPTION: The Screen-Well™ Orphan Ligand Library contains 84 compounds with defined, putative, potential or speculative activity but whose binding partners have not been identified. The library is a source of known and novel compounds for receptor de-orphaning.

CONTENTS: 84 compounds including:

D-Amino acid derivatives

Nicotine congeners

Endogenous β -carbolines

Novel Actives

Neurotransmitter metabolites

Urinary metabolites

Endo-Alkaloids

... and more

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

FORMAT: 10 mM DMSO solutions. 1 x 96-well plate. 500 μ l 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.

REFERENCES:

- 1) B. Borowsky *et al.* *Proc. Natl. Acad. Sci. USA* 2001 **98** 8966
- 2) J.R. Bunzow *et al.* *Mol. Pharmacol.* 2001 **60** 1181
- 3) A. Moser *et al.* *J. Neural Transm.* 1996 **103** 421
- 4) I.F. Musgrave *et al.* *Br. J. Pharmacol.* 2000 **129** 1057
- 5) M.M. Airaksinen *et al.* *Acta Endocrinol.* 1984 **107** 525
- 6) T.A. Green *et al.* *Biochem. Pharmacol.* 2001 **62** 1597

CITATIONS:

- 1) J.P. Saxe *et al.* *Chem. Biol.* 2007 **14** 1019

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