

## Screen-Well™ Neurotransmitter Library

**CATALOG NO.:** 2810

**Lot #:** **SAMPLE**

**DESCRIPTION:** The Screen-Well™ Neurotransmitter Library contains more than **650** CNS receptor ligands including endogenous neurotransmitters, agonists, antagonists, and marketed drugs. The library is ideal for screening or identifying recombinant orphan G protein coupled receptors, target validation, secondary screening, assay development, and other pharmacological applications.

**CONTENTS:** 650 ligands. The library contains 13 classes of ligands in 10 deep-well plates including:

- Adrenergics (Cat. # 2810)
- Cholinergics (Cat. # 2815)
- Dopaminergics (Cat. #2812)
- GABAergics (Cat. # 2819)
- Ionotropic Glutamatergics (Cat. # 2817)
- Metabotropic Glutamatergics (Cat. # 2818)
- Histaminergics and Melatonin ligands (Cat. # 2816)
- Opioids and Sigma Ligands (Cat. # 2814)
- Purinergics (Cat. # 2820)
- Serotonergics (Cat. # 2813)

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

**FORMAT:** Compounds are supplied in solution in either DMSO or water. Concentrations vary (either 10 mM or 1 mM) and are indicated for each individual compound in the product documentation. 10 x 96-well plates. 500µl per well. Each plate is also available separately.

**SUPPORTING INFORMATION:** Documentation containing a complete list of compounds and supporting information is supplied with the library or by request at [compoundlibraries@enzolifesciences.com](mailto: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.**