

SCREEN-WELL®

Neurotransmitter library (10 plate set)

The SCREEN-WELL® neurotransmitter library contains 661 CNS receptor ligands in a 96-well format. Ideal for screening or identifying recombinant orphan G protein-coupled receptors, target validation, secondary screening, validating new assays, and for routine pharmacological applications. Includes 13 classes of receptor ligands in 10 deep-well plates:

Adrenergics Histaminergics(Melatonin Ligands) Dopaminergics Ionotropic
Glutamatergics Serotonergics Metabotropic Glutamatergics Opioids(Sigma
ligands) GABAergics Cholinergics Purinergics (Adenosines)

Each ligand is supplied in solution in a biocompatible solvent at a standard concentration of 10mM (Opioid peptides are supplied at a concentrations of 1mM). Plates are available as a complete set, or individually. Supporting data are included.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-2810-0100

100 µl/well - 1Library

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	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
Contents	SCREEN-WELL® Adrenergic ligand library (BML-2811) SCREEN-WELL® Dopaminergic ligand library (BML-2812) SCREEN-WELL® Serotonergic ligand library (BML-2813) SCREEN-WELL® Opioid ligand library (BML-2814) SCREEN-WELL® Cholinergic ligand library (BML-2815) SCREEN-WELL® Histaminergic ligand library (BML-2816) SCREEN-WELL® Ionotropic Glutamatergic ligand library (BML-2817) SCREEN-WELL® Metabotropic Glutamatergic ligand library (BML-2818) SCREEN-WELL® GABAergic ligand library (BML-2819) SCREEN-WELL® Purinergic ligand library (BML-2820)
Quantity	100µl/well or 500µl/well

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com