SCREEN-WELL® Adrenergic ligand library

The adrenergic ligand library contains 83 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. The adrenergic ligand library is 1 plate of the larger neurotransmitter library (Prod. No. BML-2810) which contains over 650 CNS receptor ligands.

Citations: 3

View Online »

Ordering Information

Order Online »

BML-2811-0500	500 μl/well - 1Library
BML-2811-0100	100 μl/well - 1Library

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

one year from the date of receipt when stored at -80°C.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application HTS

Last modified: May 29, 2024



eu@enzolifesciences.com