



# SCREEN-WELL®

## Histaminergic ligand library

The Histaminergic and Melatonin Ligands Plate contains 41 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 Histaminergic and Melatonin Ligands plate is 1 plate of the larger Neurotransmitter Library (Prod. No. BML-2810) which contains over 680 ligands.

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** HTS

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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