

## A 61603 hydrobromide

A 61603 is a synthetic small molecule and potent agonist of the G protein-coupled alpha 1A receptor ( $\alpha 1A$ ), at least 35-fold selective vs  $\alpha 1B$  or  $\alpha 1D$ . More potent stimulant of phosphoinositide hydrolysis than norepinephrine, and inhibited by prazosin. Key role in modulation of the sympathetic nervous system, and increases spontaneous calcium transients more potently than phenylephrine. May be implicated in diseases such as hypertension.

### Materials Provided

|                           |                           |
|---------------------------|---------------------------|
| <b>Catalog Number:</b>    | <b>92-1230</b>            |
| <b>Quantity Supplied:</b> | 1 vial(s), 10 mg per vial |

### Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

### Product Information

|                          |   |
|--------------------------|---|
| <b>Molecular Weight:</b> | 394.79 g/mol                                    |
| <b>Source:</b>           | Synthetic                                       |
| <b>Purity:</b>           | $\geq 98\%$                                     |
| <b>Endotoxin Level:</b>  | N/A   |
| <b>Formulation:</b>      | N/A   |
| <b>Storage:</b>          | 2-8°C Please avoid multiple freeze/thaw cycles. |

### Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

|                                |                  |
|--------------------------------|------------------|
| <b>Stock Concentration:</b>    | 25 mM            |
| <b>Reconstitution Volume:</b>  | 1.013 mL         |
| <b>Reconstitution Solvent:</b> | H <sub>2</sub> O |

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)