# PNU-282987

## Nicotinic acetylcholine receptor activator

A highly selective  $\alpha 7$  nicotinic acetylcholine receptor agonist and prototypical novel therapeutic for schizophrenia.

Citations: 2

View Online »

**Ordering Information** 

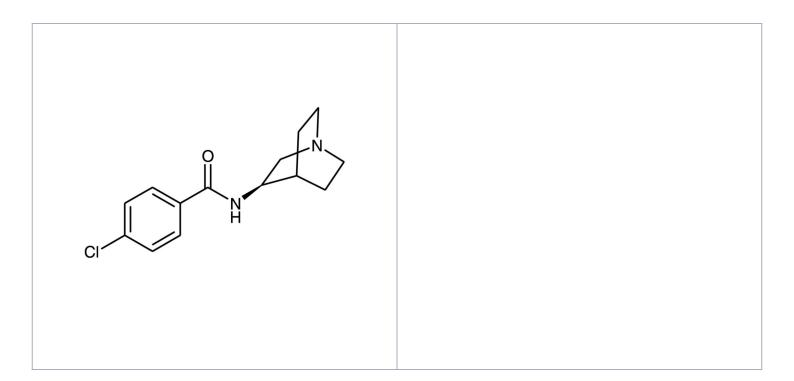
Order Online »

BML-C120-0010

10mg

Manuals, SDS & CofA

View Online »



#### **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended. Store, as supplied,

at room temperature for up to one year. Store solutions at -20°C for up to three months.

Long Term Storage Ambient

**Shipping** Ambient Temperature

### Regulatory Status RUO - Research Use Only

#### **Product Details**

Appearance White to off-white solid

**CAS** 123464-89-1

Couple Target Acetylcholine receptor

Couple Type Activator

Formula C<sub>14</sub>H<sub>17</sub>ClN<sub>2</sub>O

**MW** 264.8

Purity ≥98% (TLC)

**Solubility** Soluble in DMSO (20 mg/ml), water (6 mg/ml) or ethanol

(6 mg/ml).

Last modified: May 29, 2024