

## SC-79

AKT Activator

### Catalog # A16-901

Lot # D3370-6

CAS # 305834-79-1

### Product Description

Molecular Formula: C<sub>17</sub>H<sub>17</sub>ClN<sub>2</sub>O<sub>5</sub>

Appearance: White powder

Molecular Weight: 364.80

Purity: 98% (TLC); NMR (Conforms)

Solubilization: May be dissolved in DMSO (100 mg/ml) or ethanol (40 mg/ml, warm)

### Alias

2-Amino-6-chloro- $\alpha$ -cyano-3-(ethoxycarbonyl)-4H-1-benzopyran-4-acetic acid, ethyl ester

### Specific Activity

AKT Activator

### Storage and Stability

Store desiccated as supplied at -20°C for up to 1 year.

Store solutions at -20°C for up to 2 months.

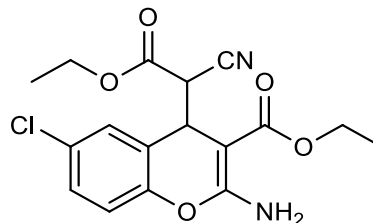
### Scientific Background

Akt activator. Binds to the pleckstrin homology domain of Akt preventing membrane translocation but activating it in the cytosol. Enhances Akt phosphorylation by upstream protein kinases. Suppresses excitotoxicity-induced neuronal death *in vitro* and *in vivo*.

### References

1. H Jo et al. *Proc. Natl. Acad. Sci. USA* 2012 109:10581

### Molecular Structure



## SC-79

AKT Activator

Catalog Number	A16-901
CAS Number	305834-79-1
Specific Lot Number	D3370-6
Purity	98% (TLC); NMR (Conforms)
Format	Lyophilized white powder
Solubilization	May be dissolved in DMSO (100 mg/ml) Ethanol (40 mg/ml, warm)
Storage & Shipping	Store desiccated as supplied at -20°C for up to 1 year. Store solutions at -20°C for up to 2 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**