

## LH-846

Casein Kinase 1 $\delta$  inhibitor

### Catalog # C67-901

Lot # D3370-5

CAS # 639052-78-1

### Product Description

Molecular Formula: C<sub>16</sub>H<sub>13</sub>ClN<sub>2</sub>OS

Appearance: White crystalline powder

Molecular Weight: 318.81

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

Solubilization: May be dissolved in DMSO (20 mg/ml) or ethanol (5 mg/ml, warm)

### Alias

N-(5-Chloro-6-methyl-2-benzothiazolyl)benzeneacetamide

### Specific Activity

Casein Kinase 1™ inhibitor

### Storage and Stability

Store desiccated as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 3 months.

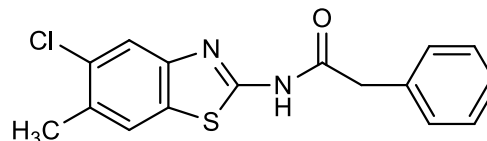
### Scientific Background

Selective inhibitor of casein kinase (CK) 1 $\delta$ , IC<sub>50</sub>s=0.29, 1.3 and 2.5  $\mu$ M for  $\delta$ ,  $\epsilon$  and  $\alpha$  isoforms respectively with no inhibition of CK2. Inhibits CK1 $\delta$ -mediated phosphorylation

### References

1. JW Lee *et al.* *Angew. Chem. Int. Ed. Engl.* 2011 50:10608

### Molecular Structure



## LH-846

Casein Kinase 1 $\delta$  inhibitor

Catalog Number	C67-901
CAS Number	639052-78-1
Specific Lot Number	D3370-5
Purity	>98% (TLC); NMR (Conforms)
Format	Lyophilized white powder
Solubilization	May be dissolved in DMSO (20 mg/ml) Ethanol (5 mg/ml, warm)
Storage & Shipping	Store desiccated as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 3 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.**