

## JNK Inhibitor IX

JNK inhibitor

### Catalog # M34-901

Lot # D3370-14

CAS # 312917-14-9

### Product Description

Molecular Formula: C<sub>20</sub>H<sub>16</sub>N<sub>2</sub>OS

Appearance: Off-White powder

Molecular Weight: 332.4

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

Solubilization: May be dissolved in DMSO (13mg/ml); or ethanol (4 mg /ml, warm)

### Alias

N-(3-cyano-4,5,6,7-tetrahydrobenzo[b]thien-2-yl)-1-naphthalenecarboxamide

### Specific Activity

JNK 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.

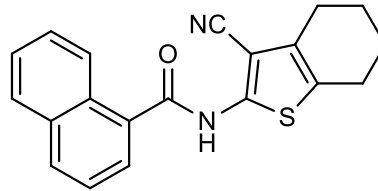
### Scientific Background

Inhibits JNK2 and JNK3 via binding to the ATP binding domain. Displays little or no activity against JNK1, p38 and a panel of 30 other kinases<sup>1</sup>.

### References

1. RM Angell *et al. Bioorg. Med. Chem. Lett.* 2007 17:1296

### Molecular Structure



## JNK Inhibitor IX

JNK inhibitor

Catalog Number M34-901  
CAS Number 312917-14-9  
Specific Lot Number D3370-14

Purity >98% (TLC); NMR (Conforms)  
Format Lyophilized powder  
Solubilization May be dissolved in DMSO (10mg/ml);  
Storage & Shipping May be dissolved in DMSO (13mg/ml); or ethanol (4 mg /ml, warm)