

## Ro 31-6045

Protein Kinase p70 (S6K) inhibitor

### Catalog # R21-901

Lot # D3370-16

CAS # 113963-68-1

### Product Description

Molecular Formula: C<sub>21</sub>H<sub>15</sub>N<sub>3</sub>O<sub>2</sub>

Appearance: Orange powder

Molecular Weight: 341.4

Purity: >98% (HPLC, TLC); NMR (conforms)

Solubilization: May be dissolved in DMSO (35 mg/ml); or ethanol (9 mg/ml)

### Alias

1-Methyl-3,4-bis(indol-3-yl)maleimide; Bisindolylmaleimide V; BIM V

### Specific Activity

Protein Kinase p70 (S6K) inhibitor

### Storage and Stability

Store desiccated as supplied at -20°C for up to 2 years.

Store solutions at -20°C for up to 3 months.

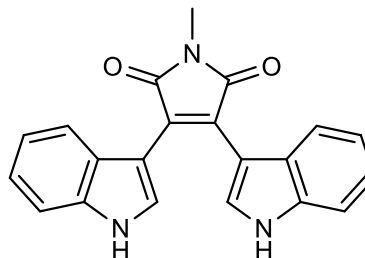
### Scientific Background

This bis indolylmaleimide derivative is a potent inhibitor of the mitogen-stimulated protein kinase p70 (S6K) acting independently of its upstream activator mTOR<sup>1</sup>. Does not inhibit PKC (at concentrations <100 μM) and therefore may also be used as a negative control for bisindolylmaleimide PKC inhibitors<sup>2</sup> such as GF 109203X<sup>3</sup> and other structurally related inhibitors<sup>4</sup>

### References

1. N Marmy-Conus et al. FEBS Lett. 2002 519:135
2. CS Harmon et al. Skin Pharmacol. 1997 10:71
3. D Toullec et al. J. Biol. Chem. 1991 266:15771
4. PD Davis et al. J. Med. Chem. 1992 35:177

### Molecular Structure



## Ro 31-6045

Protein Kinase p70 (S6K) inhibitor

Catalog #	R21-901
CAS #	113963-68-1
Lot #	D3370-16
Purity	>98% (HPLC, TLC); NMR (conforms)
Format	Orange powder
Solubilization	May be dissolved in DMSO (35 mg/ml); or ethanol (9 mg/ml)
Storage & Shipping	Store desiccated as supplied at -20°C 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.**