

## TG003

CLK family Inhibitor

### Catalog # C57-901

Lot # D3370-2

CAS # 719277-26-6

### Product Description

Molecular Formula: C<sub>13</sub>H<sub>15</sub>NO<sub>2</sub>S

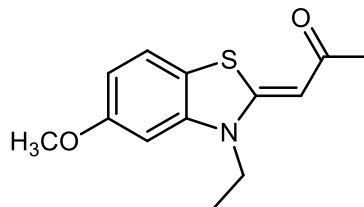
Appearance: light yellow powder

Molecular Weight: 249.33

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

Solubilization: May be dissolved in DMSO (20 mg/ml); or ethanol (15 mg/ml)

### Molecular Structure



### Alias

(Z)-1-(3-Ethyl-5-methoxy-2,3-dihydrobenzothiazol-2-ylidene)propan-2-one

### Specific Activity

CLK Family inhibitor

### Storage and Stability

Store desiccated as supplied at 4°C temperature for up to 2 years. Store solutions at -20°C for up to 2 months.

### Scientific Background

Potent, ATP-competitive inhibitor of Clk-family kinases (IC<sub>50</sub>s = 15, 20 and 200 nM for mClk4, 1 and 2 respectively and >10 μM for mClk3)<sup>1</sup>. Also inhibits DYRK1A/B (IC<sub>50</sub>s = 24 and 34 nM respectively)<sup>2</sup>.

### References

1. J Bain M Muraki *et al.* *J. Biol. Chem.* 2004 279:24246
2. A Foucourt *et al.* *Molecules* 2014 19:15411
3. S A Nishida *et al.* *Nat. Commun.* 2011 2:308

## TG003

Clk Family Inhibitor

Catalog Number	C57-901
CAS Number	719277-26-6
Specific Lot Number	D3370-2
Purity	>98% (HPLC); NMR (Conforms)
Format	Lyophilized light yellow powder
Solubilization	May be dissolved in DMSO (20 mg/ml); or ethanol (15 mg/ml)
Storage & Shipping	Store desiccated as supplied at 4°C for up to 2 years. Store solutions a -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.**