

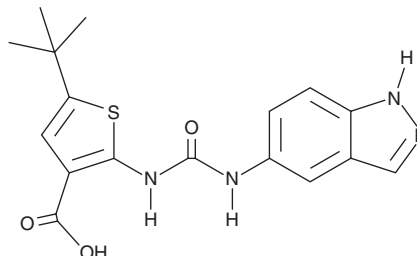
# PRODUCT INFORMATION



## S6K-18

Item No. 36624

**CAS Registry No.:** 1265789-88-5  
**Formal Name:** 5-(1,1-dimethylethyl)-2-[[[(1H-indazol-5-ylamino)carbonyl]amino]-3-thiophenecarboxylic acid  
**MF:** C<sub>17</sub>H<sub>18</sub>N<sub>4</sub>O<sub>3</sub>S  
**FW:** 358.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 225, 313 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

S6K-18 is supplied as a solid. A stock solution may be made by dissolving the S6K-18 in the solvent of choice, which should be purged with an inert gas. S6K-18 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of S6K-18 in ethanol and DMF is approximately 1 mg/ml and approximately 2 mg/ml in DMSO.

### Description

S6K-18 is an inhibitor of p70 ribosomal S6 kinase 1 (p70S6K1; IC<sub>50</sub> = 52 nM).<sup>1</sup> It is selective for p70S6K1 over a panel of 43 kinases at 1 μM.

### Reference

1. Ye, P., Kuhn, C., Juan, M., *et al.* Potent and selective thiophene urea-templated inhibitors of S6K. *Bioorg. Med. Chem. Lett.* **21**(2), 849-852 (2011).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/19/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)