

Catalogue #	Aliquot Size
<b>M16-11G -05</b>	<b>5 µg</b>
<b>M16-11G -10</b>	<b>10 µg</b>
<b>M16-11G -20</b>	<b>20 µg</b>

## COT, Active

Recombinant human protein expressed in Sf9 cells

**Catalog # M16-11G**

Lot # 1002-3

### Product Description

Recombinant human COT (30-397) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_005204](#).

### Gene Aliases

MAP3K8, EST, ESTF, TPL2, Tpl-2, c-COT, FLJ10486

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

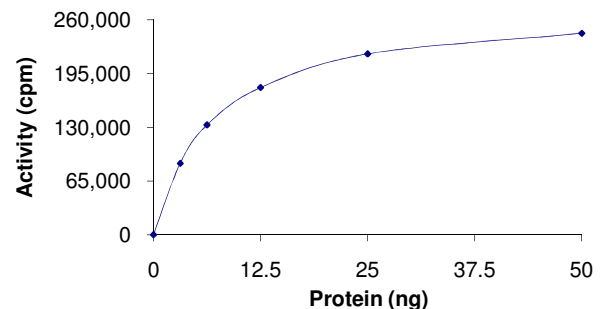
### Scientific Background

COT is an oncogene that can activate both the MAP kinase and JNK kinase pathways. COT activates I $\kappa$ B kinases and induces the nuclear production of NF- $\kappa$ B. C-terminal catalytic domain of KSR2 associates with COT and KSR2 can negatively regulate the kinase activity of COT *in vitro*. Co-transfection of KSR2 with COT in cells lead to reduced COT-mediated ERK activation and COT-induced IL8 production in a dose-dependent manner (1). COT is one of the MAP kinase kinase kinases that regulates the ERK1/ERK2 pathway in response to IL-1. Blockage of expression of COT results in failure of IL-1 to induce an increase in IL-8 and MIP-1 $\beta$  mRNA levels (2).

### References

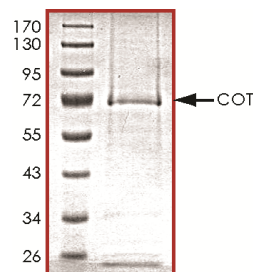
- Channavajhala, P L. et al: Identification of a novel human kinase supporter of Ras (hKSR-2) that functions as a negative regulator of Cot (Tpl2) signaling. *J. Biol. Chem.* 278: 47089-47097, 2003.
- Rodríguez C. et al: TRAF6 and Src kinase activity regulates Cot activation by IL-1. *Cell Signal.* 2006 Sep;18(9):1376-85.

### Specific Activity



The specific activity of COT was determined to be **1100 nmol /min/mg** as per activity assay protocol.

### Purity



The purity was determined to be **>75%** by densitometry. Approx. MW **~70kDa**.

## COT, Active

Recombinant human protein expressed in Sf9 cells

Catalog Number M16-11G  
Specific Activity 1100 nmol/min/mg  
Specific Lot Number I002-3

Purity >75%  
Concentration 0.1 µg/µl  
Stability 1yr At  $-70^{\circ}\text{C}$  from date of shipment  
Storage & Shipping Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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.**

# Activity Assay Protocol

## Reaction Components

### Active Kinase (Catalog #: M16-11G)

Active COT (0.1 µg/µl) diluted with Kinase Dilution Buffer III (Catalog #: K23-09) and assayed as outlined in sample activity plot. (Note: these are suggested working dilutions and it is recommended that the researcher perform a serial dilution of Active COT for optimal results).

### Kinase Dilution Buffer III (Catalog #: K23-09)

Kinase Assay Buffer I (Catalog #: K01-09) diluted at a 1:4 ratio (5X dilution) with 50ng/µl BSA solution.

### Kinase Assay Buffer I (Catalog #: K01-09)

Buffer components: 25mM MOPS, pH 7. 2, 12.5mM β-glycerol-phosphate, 25mM MgCl<sub>2</sub>, 5mM EGTA, 2mM EDTA. Add 0.25mM DTT to Kinase Assay Buffer prior to use.

### [<sup>33</sup>P]-ATP Assay Cocktail

Prepare 250µM [<sup>33</sup>P]-ATP Assay Cocktail in a designated radioactive working area by adding the following components: 150µl of 10mM ATP Stock Solution (Catalog #: A50-09), 100µl [<sup>33</sup>P]-ATP (1mCi/100µl), 5.75ml of Kinase Assay Buffer I (Catalog #: K01-09). Store 1ml aliquots at -20°C.

### 10mM ATP Stock Solution (Catalog #: A50-09)

Prepare ATP stock solution by dissolving 55mg of ATP in 10ml of Kinase Assay Buffer I (Catalog #: K01-09). Store 200µl aliquots at -20°C.

### Substrate

Unactive MEK1 (Catalog #: M02-14G) and ERK1 (Catalog #: M29-14G) were activated in a coupled reaction. Myelin Basic Protein (MBP) (Catalog #: M42-51N) diluted in distilled H<sub>2</sub>O to a final concentration of 1mg/ml was subsequently used as a substrate for the activated ERK1.

## Assay Protocol

**Step 1.** Thaw the Active COT, Kinase Assay Buffer, Unactive ERK1 and Unactive MEK1 on ice. In a pre-cooled microfuge tube, add the following reaction components bringing the initial reaction volume up to 20µl:

**Component 1.** 10µl of diluted Active COT (Catalog #M16-11G)

**Component 2.** 2µl of Unactive MEK1 (0.2µg/µl) (Catalog #M02-14G)

**Component 3.** 3µl of Unactive ERK1 (0.2µg/µl) (Catalog #M29-14G)

**Component 4.** 5µl of Kinase Dilution Buffer (Catalog #K23-09)

**Step 2.** Start the reaction by the addition of 5 µl ATP (250µM) and incubate in a water bath at 30°C for 25 minutes.

**Step 3.** After the 25 minute incubation period, remove 5µl and add to the following reaction components bringing the initial reaction volume up to 20µl:

**Component 1.** 5µl of reaction mixture

**Component 2.** 10µl distilled H<sub>2</sub>O on ice

**Component 3.** 5µl of MBP substrate (1 mg/ml) on ice (Catalog #M42-51N)

**Step 4.** Set up the blank control as outlined in step 3, excluding the addition of the substrate. Replace the substrate with an equal volume of distilled H<sub>2</sub>O.

**Step 5.** Initiate the reaction by the addition of 5µl [<sup>33</sup>P]-ATP Assay Cocktail bringing the final volume up to 25µl and incubate the mixture in a water bath at 30°C for 15 minutes.

**Step 6.** After the 15 minute incubation period, terminate the reaction by spotting 20µl of the reaction mixture onto individual pre-cut strips of phosphocellulose P81 paper.

**Step 7.** Air dry the pre-cut P81 strip and sequentially wash in a 1% phosphoric acid solution (dilute 10ml of phosphoric acid and make a 1L solution with distilled H<sub>2</sub>O) with constant gentle stirring. It is recommended that the strips be washed a total of 3 intervals for approximately 10 minutes each.

**Step 8.** Count the radioactivity on the P81 paper in the presence of scintillation fluid in a scintillation counter.

**Step 9.** Determine the corrected cpm by removing the blank control value (see Step 4) for each sample and calculate the kinase specific activity as outlined below.

### Calculation of [<sup>33</sup>P]-ATP Specific Activity (SA) (cpm/pmol)

Specific activity (SA) = cpm for 5µl [<sup>33</sup>P]-ATP / pmoles of ATP (in 5µl of a 250µM ATP stock solution, i.e., 1250 pmoles)

### Kinase Specific Activity (SA) (pmol/min/µg or nmol/min/mg)

Corrected cpm from reaction / [(SA of <sup>33</sup>P-ATP in cpm/pmol)\*(Reaction time in min)\*(Enzyme amount in µg or mg)]\*[(Reaction Volume) / (Spot Volume)]

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.**