

## ERF1 Protein

Recombinant human protein expressed in Sf9 cells

**Catalog # E380-80G**

Lot # Q2510-12

### Product Description

Recombinant human ERF1 (2-end) protein was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The ERF1 gene accession number is [NM\\_004730](#).

### Gene Aliases

eRF1; ETF1; D5S1995; ERF; RF1; SUP45L1; TB3-1

### 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°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.

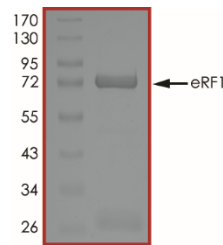
### Scientific Background

Eukaryotic peptide chain release factor subunit 1 (eRF1) directs the termination of nascent peptide synthesis (translation) in response to the termination codons UAA, UAG and UGA. Component of the transient SURF complex which recruits UPF1 to stalled ribosomes in the context of nonsense-mediated decay (NMD) of mRNAs containing premature stop codons.

### References

1. Frolova L, et al: A highly conserved eukaryotic protein family possessing properties of polypeptide chain release factor. Nature. 1994 Dec 15;372(6507):701-3.
2. <http://www.uniprot.org/uniprot/P62495>

### Purity



The purity of ERF1 was determined to be **>95%** by densitometry, approx. MW **74 kDa**.

## ERF1 Protein

Recombinant human protein expressed in Sf9 cells

|                    |   |
|--------------------|---|
| Catalog #          | E380-30G  |
| Lot #              | Q2510-12  |
| Purity             | >95%  |
| Concentration      | 0.1 µg/µl   |
| Stability          | 1yr at -70°C from date of shipment  |
| Storage & Shipping | Store product at -70°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.**

# MATERIAL SAFETY DATA SHEET

---

## Article 1 - Product Identification and Use

---

**Product Name: ERF1 Protein**

**Catalog # E380-30G**

*This product is sold only for research use by qualified laboratory personnel, and is not to be used as a drug, medical device, food additive, cosmetic, nor household chemical. It is not to be used in diagnostic, therapeutic, consumer, agricultural, nor pesticidal applications.*

Manufacturer's Name: SignalChem Pharmaceuticals Inc.  
Street Address: 110-13120 Vanier Place  
City, Prov. Postal Code: Richmond, BC, V6V 2J2  
Fax: 604-232-4601  
EMERGENCY PHONE: 604-232-4600

---

## Article 2 - Hazardous Ingredients

---

NOT AVAILABLE. We are not aware of any hazards associated with this product or its ingredients, but the chemical, physical, and toxicological properties of this product have not been investigated thoroughly. Observe normal laboratory precautions.

---

## Article 3 - Physical Data

---

This product consists of purified protein in Tris-HCl buffer shipped on dry ice. The physical properties of this product have not been investigated thoroughly.

---

## Article 4 - Fire and Explosion Hazard

---

NOT APPLICABLE

---

## Article 5 - Reactivity Data

---

NOT APPLICABLE

---

## Article 6 - Toxicologically Data

---

May be harmful by inhalation, ingestion, or skin absorption. The toxicological properties of this product have not been investigated thoroughly. Exercise due caution.

---

## Article 7 - Preventative Measures

---

Wear chemical safety goggles and compatible chemical-resistant gloves. Avoid inhalation, contact with eyes, skin or clothing.

\*\*\*\*\*MULTIPLE COMPONENT SPILL OR LEAK PROCEDURES\*\*\*\*\*

- Wear protective equipment.
  - Absorb on sand or vermiculite and place in closed containers for disposal.
  - Observe all federal, state and local environmental regulations.
- 

## Article 8 - First Aid Measures

---

- If swallowed, wash out mouth with water, provided person is conscious. Call a physician.
  - In case of skin contact, flush with copious amounts of water for at least 15 minutes. Remove contaminated clothing and shoes. If a rash or other irritation develops, call a physician.
  - If inhaled, remove to fresh air. If breathing becomes difficult, call a physician.
  - In case of eye contact, flush with copious amounts of water for at least 15 minutes while separating the eyelids with fingers. Call a physician.
- 

## Article 9 - Preparation

---

Prepared by: Jun Yan

Phone#: 1-866-954-6273

The above information is believed to be correct but does not purport to be all-inclusive and shall be used only as a guide. SignalChem shall not be held liable for any damage resulting from handling or from contact with the above product. See the Technical Specification, Packing Slip, Invoice, and Product Catalog for additional terms and conditions of sale.

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.**