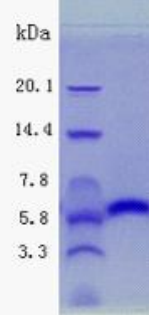


Recombinant Human EGF Protein Data Sheet

Catalog #	hRP-A0647-ES141
Size	25 µg
Protein Name	Human pro-epidermal growth factor
Protein Symbol	EGF
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	NM_001963.2
Uniprot Accession #	P01133-2
Description	This gene encodes a member of the epidermal growth factor superfamily. The encoded protein is synthesized as a large precursor molecule that is proteolytically cleaved to generate the 53-amino acid epidermal growth factor peptide. This protein acts a potent mitogenic factor that plays an important role in the growth, proliferation and differentiation of numerous cell types. This protein acts by binding the high affinity cell surface receptor, epidermal growth factor receptor. Defects in this gene are the cause of hypomagnesemia type 4. Dysregulation of this gene has been associated with the growth and progression of certain cancers. Alternate splicing results in multiple transcript variants.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	none
Peptide Length	51aa
Molecular Weight	6kDa
pI	4.3
Activity	NA
Storage	Storage buffer: 1XPBS. Store at -80°C and avoid repeated freeze-thaw cycles.
	
Reference:	• Therapeutic vaccination with an EGF-based vaccine in lung cancer: a step in the transition to a chronic disease • Epidermal growth factor (EGF) treatment on multipotential stromal cells (MSCs). Possible enhancement of therapeutic potential of MSC. • Renal magnification by EGF • Crystal structure of the complex of human epidermal growth factor and receptor extracellular domains • The EGF 61A/G polymorphism - a predictive marker for recurrence of liver metastases from colorectal cancer

GeneCopoeia™
Expressway to Discovery

GeneCopoeia Inc.

9620 Medical Center Drive, Suite 101

Rockville, MD 20850

USA

Phone: 301-762-0888

Toll free: 1-301-762-0888

Fax: 301-762-3888

Web: www.genecopoeia.com

Inquiry: inquiry@genecopoeia.com

Technical Support: support@genecopoeia.com