



DATA SHEET

Recombinant human RNase 7

Catalog no: P-400-100

Name synonyms: Ribonuclease 7, RNase7, Skin-derived antimicrobial protein 2

UniProtKB/Swiss-Prot ID Q9H1E1

Source: Human

MW: Approximately 14.5 kDa, a single non-glycosylated polypeptide chain containing 128 amino acids.

Expression host: CHO-based cell line (expressed by QMCF Technology)

Purification: Purified by affinity chromatography and gel-filtration from serum-free CHO growth media.

Concentration: 0,23 mg/ml Concentration of the protein is determined by NanoDrop using calculated Mw and extinction coefficient equation.

Buffer: 10 mM Na-acetate, pH 5.0; 150 mM NaCl

Endotoxine level: Not determined

Biological properties: Not determined

Shipping: Shipped in dry-ice.

Storage: Store at -70°C upon receipt. Recommended to aliquot into smaller quantities. Avoid repeated freeze-thaw cycles.

Related products: Monoclonal antibodies against human RNase7 For more information please visit: www.icosagen.com/antibodies

QC analysis:

Coomassie-stained SDS-PAGE

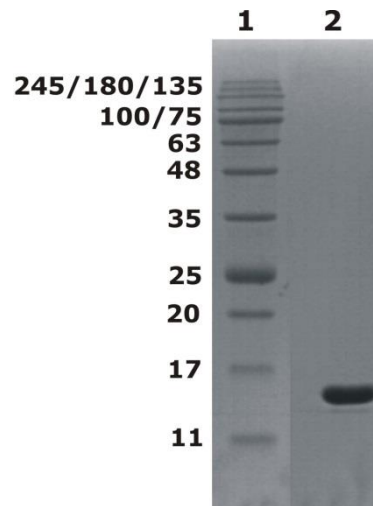


Figure 1. (A) Coomassie-stained SDS-PAGE analysis of human RNase7 Line 1. Prestained protein ladder Blue Protein Ladder Prestained, ready-to-use, 10-245 kDa (Naxo) **Line 2.** 2µg of purified RNase 7.

Background: RNase7 exhibits a potent RNase activity. Recombinant RNase 7 exhibited antimicrobial activity against uropathogens at low micromolar concentrations by disrupting the microbial membrane (Spencer et al. 2013). Has broad-spectrum antimicrobial activity against many pathogenic microorganisms and remarkably potent activity against a vancomycin resistant *Enterococcus faecium*.

This product is for research use only

distributed in the US/Canada by:
EAGLE BIOSCIENCES, INC.
20A NW Blvd, Suite 112 Nashua, NH 03063
Phone: 617-419-2019 FAX: 617-419-1110
www.EagleBio.com • info@eaglebio.com



Icosagen AS

Eerika tee 1, Õssu küla, Ülenurme vald, Tartumaa 61713, Estonia

Phone +372 737 7070, Fax +372 737 7077, info@icosagen.com

www.icosagen.com