

Lactoferrin from human milk - lyophilized viral and microbial protein

Cat. No.	Amount
PR-127	20 mg

For *in vitro* use only
Quality guaranteed for 12 months
Store at +4°C, store reconstituted aliquots at -20°C

Avoid freeze / thaw cycles

Form

Lyophilized, chromatographically purified

Source

Human milk, native

Molecular Weight

82.4 kDa

Recommended Buffer

10 mM Phosphate Buffer, 5mM NaCl

Purity

HPLC, SDS-PAGE >97%

All donors have tested negative for HIV-1, HCV antibodies and HBsAg, HIV-1 Antigen

Description

Lactoferrin is an iron binding protein. It is structurally similar to transferrin, the plasma iron transport protein. It is made by mucosal epithelium and neutrophils and is released by these cells in response to inflammatory stimuli. Bacterial growth is inhibited by its ability to sequester iron and also permeabilize bacterial cell walls by binding to lipopolysaccharides through its N-terminus. Lactoferrin can inhibit viral infection by binding tightly to the viral envelope protein. Lactoferrin appears to activate host defense systems in part by stimulating the release of interleukin-8, a neutrophil activator. It may also be involved in antibody and interleukin synthesis, lymphocyte proliferation and complement activation.

Selected References:

- B. Lonnerdaal and S.Lyer(1995) LACTOFERRIN: MOLECULAR STRUCTURE AND BIOLOGICAL FUNCTION. *AnnRev. Nutr.* **15**:93-110
- Noppe W, Plieva FM, Galaev IY, Vanhoorelbeke K, Mattiasson B, Deckmyn H. (2006) Immobilised peptide displaying phages as affinity ligands. Purification of lactoferrin from defatted milk. *J. Chromatogr A.* **6**:1101(1-2):79-85
- Legrand D, Elass E, Carpentier M, Mazurier (2006) Interactions of lactoferrin with cells involved in immune function. *J. Biochem Cell Biol.* **84**(3):282-90
- Seganti L, Di Biase AM, Marchetti M, Pietrantonio A, Tinari A, Superti F. (2004) Antiviral activity of lactoferrin towards naked viruses. *Biomaterials.* **17**(3):295-9.
- van der Strate BW, Beljaars L, Molema G, Harmsen MC, Meijer DK. (2001) Antiviral activities of lactoferrin. *Antiviral Res.* **52**(3):225-39.