

---

## DMKN Polyclonal Antibody, FITC Conjugated

(Catalog # A58812)

---

### Background

May act as a soluble regulator of keratinocyte differentiation.

### Description

DMKN Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q6E0U4

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Dermokine protein (1-137AA)

### Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Dermokine, DMKN, UNQ729/PRO1411, Epidermis-specific secreted protein SK30/SK89

### Application

ELISA; Recommended dilution: ELISA 1:100-1:500

---