

# PRODUCT INFORMATION

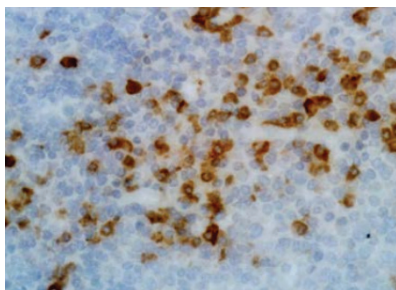


## Igk Light Chain (human) Rabbit Monoclonal Antibody (Clone RM126) Item No. 32108

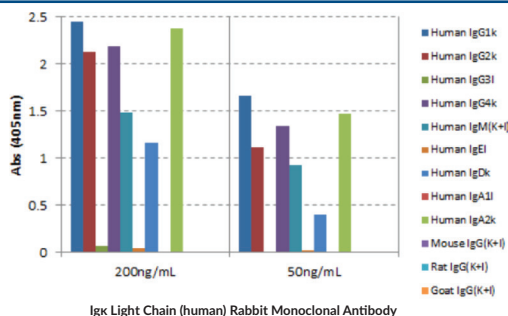
### Overview and Properties

**Contents:** This vial contains 100 µg of protein A-affinity purified monoclonal antibody.  
**Synonym:** Immunoglobulin κ Light Chain  
**Immunogen:** Human IgG  
**Cross Reactivity:** (+) Igk light chain; (-) Human λ light chain, cynomolgus monkey, goat, mouse, rat, rhesus monkey IgG  
**Species Reactivity:** (+) Human  
**Form:** Liquid  
**Storage:** -20°C (as supplied)  
**Stability:** ≥1 year  
**Storage Buffer:** PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide  
**Concentration:** 1 mg/ml  
**Clone:** RM126  
**Host:** Rabbit  
**Isotype:** IgG  
**Applications:** ELISA, flow cytometry (FC), immunocytochemistry (ICC), and immunohistochemistry (IHC); the recommended starting concentration for ELISA is 0.05-0.2 µg/ml and 0.5-2 µg/ml for FC, ICC, and IHC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

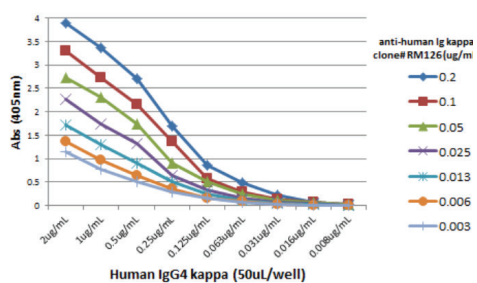
### Images



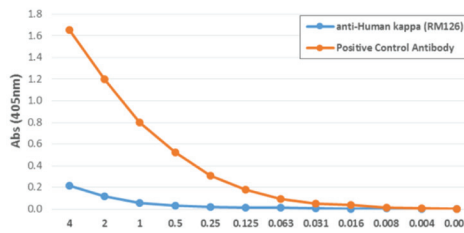
Immunohistochemical staining of formalin-fixed and paraffin-embedded human tonsil tissue section using Igk Light Chain (human) Rabbit Monoclonal Antibody (Clone RM126).



ELISA of Human Immunoglobulins (Igs). Igk Light Chain (human) Rabbit Monoclonal Antibody (Clone RM126) reacts only to the κ light chain of human Igs and not to λ light chain, mouse IgG, rat IgG, or goat IgG.



A Titer ELISA using Igk Light Chain (human) Rabbit Monoclonal Antibody (Clone RM126). The plate was coated with different amounts of human IgG4κ. A serial dilution of Igk Light Chain (human) Rabbit Monoclonal Antibody (Clone RM126) was used as the primary antibody, followed by an alkaline phosphatase conjugated anti-rabbit IgG as the secondary antibody.



ELISA of Rhesus Monkey IgG. Igk Light Chain (human) Rabbit Monoclonal Antibody (Clone RM126) does not react to monkey IgG. The plate was coated with rhesus monkey IgG. A serial dilution of Igk Light Chain (human) Rabbit Monoclonal Antibody (Clone RM126) and a monkey IgG binding antibody (positive control) was used as the detection antibody.

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/09/2023

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

Ig $\kappa$  light chain is one type of light chain found in immunoglobulins, which are part of the immunoglobulin superfamily of glycoproteins that plays a central role in the adaptive immune response.<sup>1</sup> Immunoglobulins are produced by B cells and later secreted by plasma cells as antibodies.<sup>2</sup> They are composed of two heavy chains of approximately 50 kDa each and two light chains of approximately 25 kDa each.<sup>1</sup> The heavy chains are linked together by disulfide bonds to form an Fc region and also combine with the light chains to form the Fab region, which mediate receptor and antigen binding, respectively.<sup>3</sup> Mammalian immunoglobulins contain either Ig $\kappa$  or Ig $\lambda$  light chains, each of which are composed of a constant and variable domain.<sup>4</sup> The ratio of Ig $\kappa$  to Ig $\lambda$  light chain-containing antibodies varies between species, with ratios of 20:1, 2:1, and 1:20 in mice, humans, and cattle, respectively. Ig $\kappa$  and Ig $\lambda$  free light chains (FLCs) are produced during immunoglobulin synthesis, and accumulation of these FLCs, primarily Ig $\kappa$ , is associated with various disorders, including light-chain deposition disease, multiple myeloma, rheumatoid arthritis, diabetic nephropathy, and systemic lupus erythematosus (SLE).<sup>2,5,6</sup> Cayman's Ig $\kappa$  Light Chain (human) Rabbit Monoclonal Antibody (Clone RM126) can be used for ELISA, flow cytometry (FC), immunocytochemistry (ICC), and immunohistochemistry (IHC) applications. The antibody recognizes the  $\kappa$  light chain from human samples.

## References

---

1. Schroeder, H.W., Jr. and Cavicini, L. Structure and function of immunoglobulins. *J. Allergy Clin. Immunol.* **125**(2 Suppl. 2), S41-S52 (2010).
2. Esparvarinha, M., Nickho, H., Mohammadi, H., *et al.* The role of free kappa and lambda light chains in the pathogenesis and treatment of inflammatory diseases. *Biomed. Pharmacother.* **91**, 632-644 (2017).
3. Vaillant A.A.J. and Ramphul K. Immunoglobulin. In: StatPearls [Internet]. Treasure Island (FL): StatPearls Publishing (2020). Available from: <https://www.ncbi.nlm.nih.gov/books/NBK513460/>
4. Janeway, C.A., Jr., Travers, P., Walport, M., *et al.* Antigen recognition by B-cell and T-cell receptors. *Immunobiology*. 6th edition, *Garland Science Publishing* (2004).
5. Jimenez-Zepeda, V.H. Light chain deposition disease: Novel biological insights and treatment advances. *Int. J. Lab. Hematol.* **34**(4), 347-355 (2012).
6. Sannier, A., Hanouna, G., Daugas, E., *et al.* IgA kappa light and heavy chain deposition disease in multiple myeloma. *Br. J. Haematol.* **183**(1), 13 (2018).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM