

# Protein L Biosensors

For whole molecule antibody and antibody fragment quantitation

## Key features

- Broader recognition of Ig isotypes and species than Protein A or G
- Regenerable and cost-effective format
- Binds whole molecule Ab and FAb fragments
- Quantitate in crude matrices

## Overview

Dip and Read™ Protein L (ProL) biosensors provide a rapid and direct method for quantifying a broad set of kappa light chain containing immunoglobulins, including whole molecules, FAb fragments and single chain variable fragments, in buffer, conditioned media or complex matrices. Protein L, which is factory-immobilized onto the biosensor, binds antibodies through the kappa light chain and recognizes a wider range of antibody classes than either Protein A or Protein G, including IgG, IgM, IgA, IgE and IgD. Due to its specificity, Protein L recognizes mouse Vkl (but not VkII, VkIII and VkIV) and human Vkl, VkIII and VkIV (but not VkII). The Protein L Biosensor is especially useful for quantifying antibodies and antibody fragments from serum based cultures because Protein L does not bind bovine immunoglobulins, which often contaminate serum supplements.

## Quick facts

- **Typical dynamic range for hFAB:**
  - 0.05–2000 µg/mL (Octet® RED96 and RED384 systems)
  - 0.05–700 µg/mL (Octet QK® and QK384 systems)
  - 0.1–700 µg/mL (Octet QK system)
- Sensitivity can be extended by longer read times and higher shake speed.
- **Throughput:**
  - Up to 384 samples < 60 minutes (Octet RED384 system)
  - Up to 96 samples < 30 minutes (Octet RED96 system)
- **Precision:** ≤ 10% CVs

## Principle of antibody quantitation

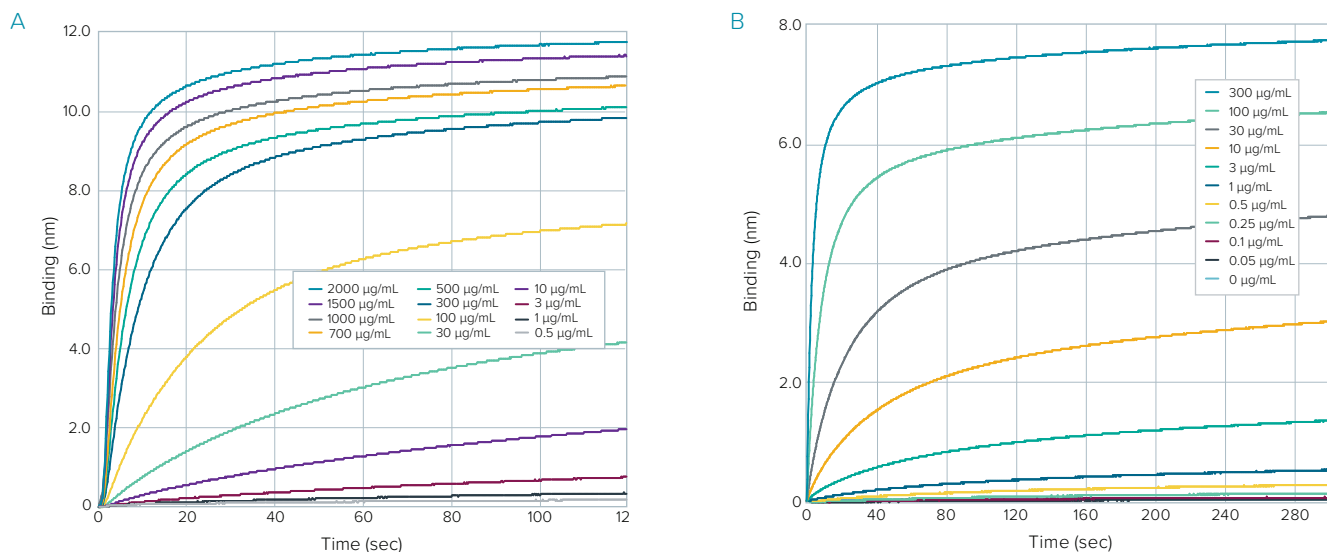
Protein L is an immunoglobulin binding protein originally isolated from *Peptostreptococcus magnus* that binds the kappa light chains of most antibodies and antibody fragments. Immunoglobulin

binding to the Protein L Biosensor alters the properties of light on the biosensor surface, allowing the association event to be monitored in real time using the Octet system. Higher antibody concentrations cause faster binding rates. Analyte concentration is extrapolated from a standard curve. Due to the high specificity of Protein L for the antibody light chain, Protein L biosensors can directly quantitate analytes in conditioned media and complex matrices.

Antibody	Protein L	Protein A	Protein G
Rat Total IgG	s	w	m
Rat IgG1	s	w	m
Rat IgG2a	s	nb	s
Rat IgG2b	s	nb	w
Rat IgG2c	s	s	s
Rabbit IgG	w	s	s
Mouse IgG1	s	w	m
Mouse IgG2a	s	s	s
Mouse IgG2b	s	s	s
Mouse IgG3	s	s	s
Goat IgG	nb	w	s
Sheep IgG	nb	w	s
Bovine IgG	nb	w	s
Guinea Pig IgG	w	s	w
Hamster IgG	s	m	m
Pig IgG	s	s	w
Horse IgG	–	w	s
Donkey IgG	–	m	s
Dog IgG	–	s	w
Cat IgG	–	s	w
Human IgG1	s	s	s
Human IgG2	s	s	s
Human IgG3	s	w	s
Human IgG4	s	s	s

s = strong m = medium w = weak nb = no binding

**Table 1:** Proteins A, G and L possess unique affinity profiles toward antibodies from different species and isotypes. Protein L binds with high affinity to several more species and isotypes than Proteins A and G, broadening its potential applications.



**Figure 1:** Protein L detects a human IgG FAb over a wide dynamic range from 0.5–2000 µg/mL in sample diluent using parameters tuned for high (A) and low (B) concentrations. High concentration samples (0.5–2000 µg/mL, Panel A) were shaken at 400 rpm. Low concentration samples (0.05–300 µg/mL, Panel B) were shaken at 1000 rpm.

Biosensor #	Concentration (µg/mL) of hFAB detected per regeneration cycle										Statistics			
	R0	R1	R2	R3	R4	R5	R6	R7	R8	R9	Avg.	Std. Dev.	CV%	% Loss
1	10.2	10.1	10.0	10.0	10.0	10.0	9.91	9.87	9.92	9.85	9.99	0.11	1.1%	-3.4%
2	10.3	10.2	10.1	10.1	10.1	10.1	10.0	9.96	10.0	9.90	10.1	0.12	1.2%	-3.9%
3	10.4	10.3	10.3	10.3	10.2	10.2	10.2	10.1	10.1	10.1	10.2	0.10	1.0%	-2.9%

**Table 2:** Three Protein L biosensors were used to repeatedly analyze samples of 10 µg/mL human IgG FAb in neat CHO DG44 medium at a shake rate of 400 rpm. The Protein L biosensors were pre-conditioned with 3 regeneration cycles prior to the first measurement (R0) and regenerated in 10 mM Glycine pH 1.5 between each subsequent measurement (R0 through R9). After 10 regeneration cycles, the loss of capacity was less than 4% and the coefficient of variance was less than 2%, demonstrating efficient recovery of capacity and high reproducibility.

## Wide dynamic range

For most human antibodies and antibody fragments, Protein L biosensors detect analytes at concentrations from 50 ng/mL to 2 mg/mL with variation of the assay read time and shake speed (Figure 1). The ability to tune the assay over this wide dynamic range reduces the need for multiple dilutions, increasing the number of samples per plate and assay efficiencies.

## Cost-effective regeneration

Protein L biosensors can be regenerated up to ten times or more via a standard low-pH protocol. Regeneration dissociates the bound antibody from the Protein L Biosensor, allowing additional analyses and providing a cost-effective format for analyzing large sample libraries. Table 2 summarizes the high recovery and precision of three representative biosensors over ten regeneration cycles.

## Ordering information

Part no.	UOM	Description
18-5085	Tray	96 biosensors coated with Protein L (recommended Sample Diluent sold separately)
18-5086	Pack	Five trays of 96 biosensors coated with Protein L (recommended Sample Diluent sold separately)
18-5087	Case	Twenty trays of 96 biosensors coated with Protein L (recommended Sample Diluent sold separately)

**Note:** A purified standard that is identical to the experimental samples is required.

Protein A and Protein G biosensors are also available.



**ForteBio**  
47661 Fremont Boulevard  
Fremont, CA 94538  
888.OCTET-75 or 650.322.1360  
fortebio.info@moldev.com

**ForteBio Analytics (Shanghai) Co., Ltd.**  
No. 88 Shang Ke Road  
Zhangjiang Hi-tech Park  
Shanghai, China 201210  
salesops.china@moldev.com

**Molecular Devices (UK) Ltd.**  
660-665 Eskdale  
Winnersh Triangle  
Wokingham, Berkshire  
RG41 5TS, United Kingdom  
+44 118 944 8000  
uk@moldev.com

**Molecular Devices (Germany) GmbH**  
Bismarckring 39  
88400 Biberach an der Riss  
Germany  
+00800 665 32860