

1 Image

Product name	Feline IL-8 ELISA Kit
---------------------	-----------------------

Detection method	Colorimetric
------------------	--------------

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

Sample type	Cell culture supernatant, Serum, Plasma
--------------------	---

Assay type	Sandwich (quantitative)
-------------------	-------------------------

Sensitivity 25 pg/ml

Range	25 pg/ml - 6000 pg/ml
--------------	-----------------------

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	89.71	82% - 97%
Plasma	84.6	71% - 95%
Cell culture media	118	100% - 149%

Assay duration	Multiple steps standard assay
-----------------------	-------------------------------

Species reactivity **Reacts with:** Cat

Product overview Abcam's IL-8 Feline ELISA Kit (ab193688) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of feline IL-8 in serum, plasma (feline IL8 concentration is pretty low in normal plasma, it may not be detected in this assay) and cell culture supernatant.

This assay employs an antibody specific for feline IL-8 coated on a 96-well plate. Standards and

samples are pipetted into the wells and the immobilized antibody captures IL-8 present in the samples. The wells are washed and biotinylated anti-feline IL8 antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of IL-8 bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

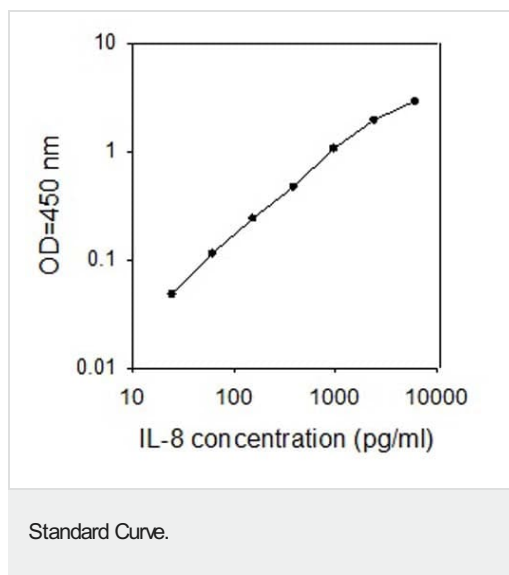
Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Biotinylated feline IL-8 detection antibody	2 vials
feline IL-8 Standards (lyophilized)	2 vials
Pre-coated feline IL-8 Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Relevance IL-8 is a chemotactic factor that attracts neutrophils, basophils, and T-cells, but not monocytes. It is also involved in neutrophil activation. It is released from several cell types in response to an inflammatory stimulus. IL-8(6-77) has a 5-10-fold higher activity on neutrophil activation, IL-8(5-77) has increased activity on neutrophil activation and IL-8(7-77) has a higher affinity to receptors CXCR1 and CXCR2 as compared to IL-8(1-77), respectively.

Cellular localization Secreted

Images



Standard Curve using ab193688 IL-8 Feline ELISA Kit.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors