

Nacka Strand, September 1, 2016



## Launch of new Canine IFN- $\gamma$ ELISA and ELISpot kits

Dear Researcher,

We are pleased to launch new kits and reagents for detection of canine IFN- $\gamma$  as of September 29, 2016.

Mabtech has developed novel monoclonal antibodies (mAbs) reactive with canine IFN- $\gamma$ . These mAbs will be included in our ELISA and ELISpot kits and available as separate items. The antibodies currently included in these kits will be discontinued. The performance of the new mAbs is highly comparable with the discontinued. Please see the back of this page for a list of all our canine IFN- $\gamma$  products.

If you have any questions regarding our products we encourage you to contact us and we welcome any comments you might have.

A handwritten signature in black ink, appearing to read "Niklas Ahlborg".

Niklas Ahlborg, Assoc. Prof.  
Director of Research and Development, Mabtech

[www.mabtech.com](http://www.mabtech.com)

### Contact Details

Email:	<a href="mailto:mabtech.usa@mabtech.com">mabtech.usa@mabtech.com</a>	Mabtech, Inc.
Phone:	+1 513 871 4500	M.E.B Suite 220, 3814 West Street
Fax:	+1 513 871 7353	OH 45227 Cincinnati
		United States

## Monoclonal antibodies

anti-canine IFN- $\gamma$ mAb MT13, purified	250 $\mu$ g	3113-3-250
	1 mg	3113-3-1000
anti-canine IFN- $\gamma$ mAb MT166, biotinylated	250 $\mu$ g	3113-6-250
	1 mg	3113-6-1000

## ELISA

Canine IFN- $\gamma$ ELISA development kit (ALP)	for 6 plates	3113-1A-6
	for 20 plates	3113-1A-20
Canine IFN- $\gamma$ ELISA development kit (HRP)	for 6 plates	3113-1H-6
	for 20 plates	3113-1H-20

Ready-to-use pNPP and TMB ELISA substrates are available in our web shop.

## ELISpot

Canine IFN- $\gamma$ ELISpot <sup>BASIC</sup> (ALP)	for 4 plates	3113-2A
Canine IFN- $\gamma$ ELISpot <sup>BASIC</sup> (HRP)	for 4 plates	3113-2H
Canine IFN- $\gamma$ ELISpot <sup>PLUS</sup> (ALP)	2 plates	3113-4APW-2
	10 plates	3113-4APW-10
Canine IFN- $\gamma$ ELISpot <sup>PLUS</sup> (HRP)	2 plates	3113-4HPW-2
	10 plates	3113-4HPW-10

Please note that the product codes for the ELISA and ELISpot products are the same as before.