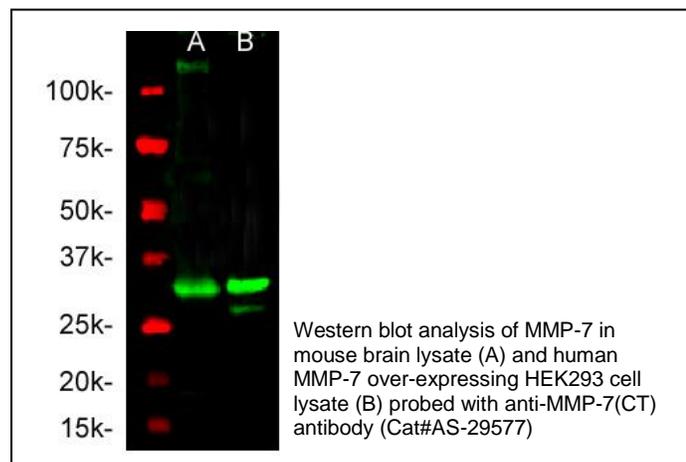


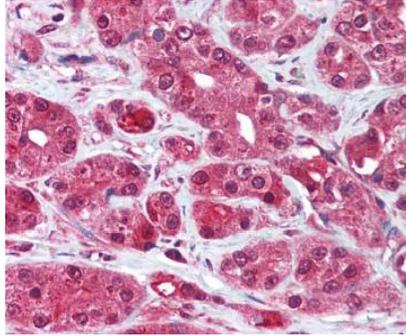
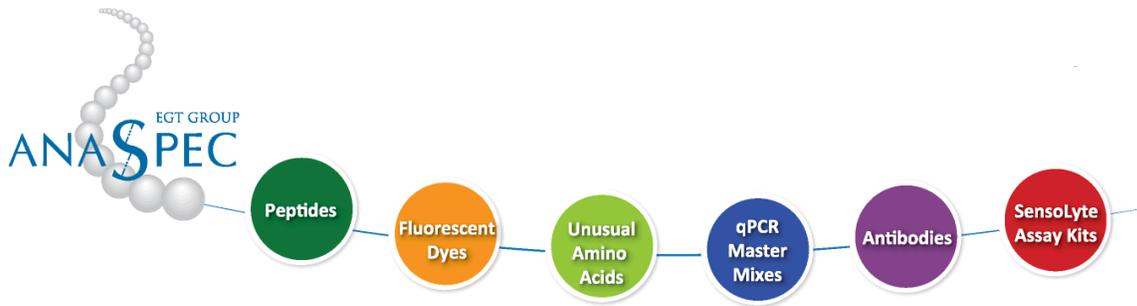


Update: June, 2017

## Product Data Sheet

<b>Product Name:</b>	Anti-MMP-7 (CT) <i>Rabbit Polyclonal Antibody</i>
<b>Catalog Number:</b>	AS-29577
<b>Lot Number:</b>	See label on vial
<b>Storage Buffer:</b>	1X PBS (pH 7.4) containing 0.05% sodium azide and <0.1% BSA
<b>Size:</b>	50 µg
<b>Concentration:</b>	0.2 mg/mL
<b>Immunogen:</b>	KLH conjugated with synthetic peptide corresponding to the C-terminus of human MMP-7.
<b>Species Reactivity:</b>	Species reactivity includes human and mouse while others remain unknown.
<b>Application Notes:</b>	The following concentration ranges are recommended starting points for this product. Optimal working concentrations should be determined by the investigator for specific applications.
	Western blot: 0.5-2.0 µg/mL
	Immunohistochemistry: 10.0-20.0 µg/mL





Immunocytochemistry analysis of FFPE human carcinoma tissue stained with anti-MMP-7(CT) antibody (Cat#AS-29577).

**Background:**

Matrix metalloproteinases (MMPs) belong to a family of proteases that are essential for the breakdown of extracellular matrix (ECM). They are thus important in apoptosis; tumor cell growth, invasion, and metastasis; as well as angiogenesis and wound healing.<sup>1-3</sup> Most MMPs contain common domain structures that include a signal sequence, a propeptide, a catalytic domain, and a hemopexin-like (Hpx) domain.<sup>4</sup> MMP-7, also known as matrilysin, is responsible for hydrolyzing proteoglycans and ECM glycoproteins.<sup>4</sup> It is expressed in normal cells, specifically glandular epithelial cells.<sup>5</sup>

**Storage:**

Store at 4 °C for 1-2 weeks. Aliquot and store at -20°C up to 1 year. Avoid freeze and thaw cycle.

**References:**

1. Itoh, Y. et al. *Essays in Biochemistry* **38**, 21 (2002).
2. Ayake, T. et al. *Nature Immunol* **1**, 113 (2000).
3. Ito, A. et al. *J Biol Chem* **271**, 14651 (1996).
4. Fowlkes, J. et al. *J Biol Chem* **270**, 274 (1995).
5. Woessner, J. et al. *FASEBJ* **5**, 2145 (1991).

This product is for *in vitro* research use only.