

## TFF1 (trefoil factor 1 or pS2 protein) Monoclonal Antibody

Cat No: G01094 - 100  $\mu$ L

### General Data

<b>Shipping:</b>	wet ice
<b>Stability:</b>	Store at -20degrees; shelf life 3 year maximum after production Avoid repeat freezing and thawing cycles. Does not contain any preservative
<b>Formulation:</b>	Liquid Ascite, does not contain any preservative therefore avoid repeat freezing and thawing cycles
<b>Host:</b>	Mouse
<b>Antigen:</b>	Synthetic peptide
<b>Clone:</b>	3D5
<b>Isotype:</b>	IgG1, $\kappa$



<b>Application(s):</b>	ELISA
	Western Blot
	Immunocytochemistry
	Immunohistochemistry
	Recommended dilutions: 1/1000-1/10000

**Specificity:** human, does not cross react on mouse

### Scientific Literature

Rio MC, Bellocq JP, Daniel JY, Tomasetto C, Lathe R, Chenard MP, Batzenschlager A, Chambon P. Breast cancer-associated pS2 protein: synthesis and secretion by normal stomach mucosa. Science 1988; 241: 705-708.

FP/10/20

For research laboratory use only – Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

#### Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac  
10 bis avenue Ampère  
78180 Montigny le Bretonneux - France  
Tel.: +33 (0)139 306 036

<https://www.bertin-bioreagent.com/pa206/contact-us>  
<https://www.bertin-bioreagent.com>