

# Trefoil factor 2 / TFF2 Monoclonal Mouse Antibody (GE16C)



## Product Description

Trefoil peptides are protease resistant molecules secreted throughout the gut that play a role in mucosal healing. These peptides contain three intrachain disulfide bonds, forming the trefoil motif, or P-domain.

This antibody is available purified, with BSA and azide (0.2 mg/mL) or purified, BSA- and azide-free (1 mg/mL).

## References

Note: References for this clone sold by other suppliers may be listed for expected applications.

Clin Cancer Res (2005) 11(18): 6472-6478. (IHC, FFPE)

### Product attributes

Antibody number	#3060
Antibody reactivity (target)	TFF2, Trefoil factor 2
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	GE16C
Isotype	IgM, kappa
Molecular weight	14 kDa
Synonyms	SML1; SP; Spasmolysin; Spasmodytic polypeptide; TFF2; trefoil factor 2 (spasmodytic protein 1); Trefoil factor 2; Trefoil factor 2 precursor
Human gene symbol	TFF2
Entrez gene ID	7032
SwissProt	Q03403
Unigene	2979
Immunogen	Synthetic peptide corresponding to a 16-amino acid C-terminal sequence of the human Trefoil factor 2.
Antibody target cellular	Secreted (extracellular)
Species reactivity	Human
Expected antibody	IHC (FFPE) (published for clone)
Antibody application notes	ELISA: For coating, order Ab without BSA & azide; Optimal dilution for a specific application should be determined. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
Positive control	Human stomach or duodenum tissues (IHC).
Shipping condition	Room temperature
Storage Conditions	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended

This datasheet was generated on September 2, 2024 at 12:24:50 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/trefoil-factor-2-tff2-monoclonal-mouse-antibody-ge16c/>