

# GCDFP-15 (human) monoclonal antibody (D6)

Gross cystic disease is a common premenopausal disorder in which gross cysts are the predominant pathologic lesion. It is characterized by production of a fluid secretion which accumulates in the breast cysts. Gross cystic disease fluid is a pathologic secretion from breast composed of several glycoproteins, including a unique 15kDa monomer protein, GCDFP-15. The cells within the body that produce GCDFP-15 appear to be restricted primarily to those with apocrine function. Studies have found GCDFP-15 to be a highly specific and sensitive marker for breast cancer. GCDFP-15 is expressed by apocrine sweat glands, eccrine glands (variable), minor salivary glands, bronchial glands, metaplastic epithelium of the breast, benign sweat gland tumors of the skin and the serous cells of the submandibular gland. Breast carcinomas (primary and metastatic lesions) with apocrine features express GCDFP-15. GCDFP-15 is positive in extramammary Paget's Disease. Other tumors tested are negative.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	BRST-2
Application	IHC (PS)
Application Notes	This antibody is heat sensitive. Paraffin sections that have been heat fixed at 60°C may lose antibody binding site.
Clone	D6
Formulation	Liquid. In 0.01M PBS containing 1% BSA and 0.1% sodium azide .
Host	Mouse
Immunogen	GCDFP-15 (gross cystic disease fluid protein-15; BRST-2).
Isotype	IgG2a
Positive Control	Human axilla or breast carcinoma.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections): Formalin-fixed, paraffin-embedded tissues; pre-treatment not required; suggested dilution: ≥1:40 for biotin based detection. For optimal staining, incubate primary antibody 20-60 minutes at room temperature). Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P12273

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)