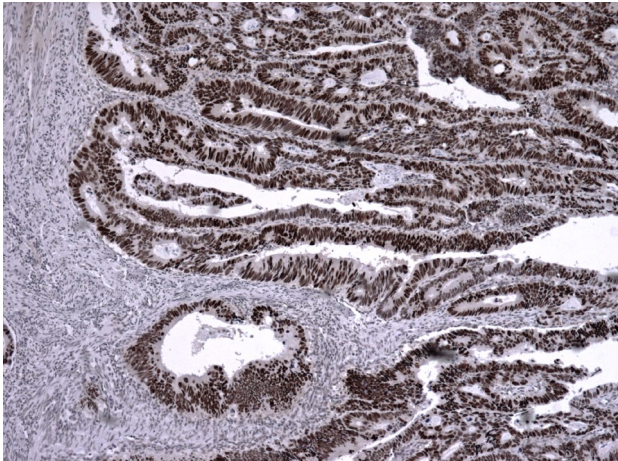


Mouse monoclonal anti-p53 antibody, clone BP-53-12

For Research Use Only

Cat. №	Volume	Configuration	Approx. number of tests
10-310009-7	7 ml	Pre-diluted Antibody, Ready-To-Use	50
10-310009-01	0.1 ml	Concentrated Antibody	75
10-310009-05	0.5 ml		375
10-310009-1	1 ml		750

Intended Use



RUO — Research Use Only.

This product is used to qualitatively detect p53 in normal and neoplastic formalin fixed paraffin embedded tissue sections in immunohistochemical detection methodology.

Description

Antibodies bind to the normal and mutant types of p53 protein and are used to detect increased expression of p53. For many types of tumors, the detection of p53 is predictive. The

wild-type p53 protein is present in normal cells, but, due to the extremely short half-life, it is not detected using standard immunohistochemical methods. Somatic mutations of the p53 gene are often found in various tumor processes. The mutant p53 protein is more stable than the wild-type protein and accumulates in tumor cells in significant concentration.

Antigen

N-terminal part of recombinant protein

Reagent provided

This antibody is purified immunoglobulin IgG2a diluted in 10 mM Phosphate buffered saline (PBS), pH 7.4 containing 1% bovine serum albumin (BSA) and 0.05% ProClin™ 300 as antimicrobial agent.

Uses/Limitations:

For professional users.

For Research Use Only.

Do not use if reagent becomes cloudy.

Do not use past expiration date.



Contains ProClin™ 300 as a preservative (0.05% w/v)

Usage dilution

10-310009-7: Ready to Use

10-310009-01; 10-310009-05; 10-310009-1: Dilute 1:50 to 1:100 with Antibody Diluent (REF 12-310001) before use. Optimum dilution factor may vary depending on the specimen and preparation process and should be determined by each individual investigator.

Epitope retrieval

Staining of formalin fixed, paraffin embedded tissue sections is significantly enhanced by pretreatment with Citrate Buffer, pH 6.0 (REF 12-310002)

Staining procedure

Incubate this antibody with tissue section for 30-60 minutes at room temperature. Follow the instructions from the selected detection system.

Positive control tissue

Colon adenocarcinoma

Staining pattern

Nucleus

Storage

Store at 2-8°C and in the dark. Do not use after expiration date.

Warning and precautions



Refer to SDS.

Support

Please contact ClearMe LLC by phone or by e-mail.

info@diagomix.com

Symbols

REF	LOT		
Catalog number	Batch number	Temperature Range	Use by

