



Catalog No. 09-0002

## HER2 Control Slides

### EXPECTED STAINING:

Lot No.

MARKER	IMMUNOSTAIN PATTERN
HER2 (c-erbB-2)	Membrane

Positive tissue controls should only be utilized for monitoring the correct performance of processed tissues and test reagents, rather than as an aid in the interpretation of samples. If the positive tissue controls fail to demonstrate positive staining, results with the test specimens should be considered invalid.

### TISSUE PROVIDED:

Source: Breast carcinoma  
Fixative: 10% Neutral buffered formalin  
Embedding: Paraffin  
Thickness: 4-6 µm  
Mounting: Mounted on positively charged slides and baked for 15 min. at 59°C.

**QUANTITY:** 5 unstained slides.

**STORAGE:** 2-8°C.

**INTENDED USE:** For research use only. Deparaffinize and rehydrate before use. Process positive control slides using same protocol as test samples.

### LIMITATIONS:

- Immunohistochemistry is a multistep diagnostic process that consists of specialized training in the selection of the appropriate reagents; tissue selection, fixation, and processing; preparation of the IHC slide; and interpretation of the staining results. It is the responsibility of a qualified individual who is familiar with the proper use of IHC antibodies, reagents and methods to interpret all of the steps used to prepare and interpret the final IHC preparation.
- The possibility of unexpected reactions even in tested tissue cannot be completely eliminated due to biological variability of antigen expression in normal and abnormal tissues. Contact Invitrogen at 800.955.6288 with documented unexpected reaction(s).
- False-positive results may be seen due to non-immunological binding of proteins or substrate reaction products. They may also be caused by pseudoperoxidase activity (erythrocytes), endogenous peroxidase activity (cytochrome c), or endogenous biotin (e.g. liver, breast, brain, kidney) (depending on the type of immunostain used).
- This tissue section is not known to contain any infectious agents, however, standard precautions should be observed.

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288

E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PIN 30693  
1089

(Rev 08/08) DCC-08-

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.



Catalog No. 09-0002

## HER2 Control Slides

### EXPECTED STAINING:

Lot No.

MARKER	IMMUNOSTAIN PATTERN
HER2 (c-erbB-2)	Membrane

Positive tissue controls should only be utilized for monitoring the correct performance of processed tissues and test reagents, rather than as an aid in the interpretation of samples. If the positive tissue controls fail to demonstrate positive staining, results with the test specimens should be considered invalid.

### TISSUE PROVIDED:

Source: Breast carcinoma  
Fixative: 10% Neutral buffered formalin  
Embedding: Paraffin  
Thickness: 4-6 µm  
Mounting: Mounted on positively charged slides and baked for 15 min. at 59°C.

**QUANTITY:** 5 unstained slides.

**STORAGE:** 2-8°C.

**INTENDED USE:** For research use only. Deparaffinize and rehydrate before use. Process positive control slides using same protocol as test samples.

### LIMITATIONS:

- Immunohistochemistry is a multistep diagnostic process that consists of specialized training in the selection of the appropriate reagents; tissue selection, fixation, and processing; preparation of the IHC slide; and interpretation of the staining results. It is the responsibility of a qualified individual who is familiar with the proper use of IHC antibodies, reagents and methods to interpret all of the steps used to prepare and interpret the final IHC preparation.
- The possibility of unexpected reactions even in tested tissue cannot be completely eliminated due to biological variability of antigen expression in normal and abnormal tissues. Contact Invitrogen at 800.955.6288 with documented unexpected reaction(s).
- False-positive results may be seen due to non-immunological binding of proteins or substrate reaction products. They may also be caused by pseudoperoxidase activity (erythrocytes), endogenous peroxidase activity (cytochrome c), or endogenous biotin (e.g. liver, breast, brain, kidney) (depending on the type of immunostain used).
- This tissue section is not known to contain any infectious agents, however, standard precautions should be observed.

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288

E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PIN 30693  
1089

(Rev 08/08) DCC-08-

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.