

p185(HER-2) antibody, FITC (mAb)

Mouse Anti-Human p185 (Human EGF Receptor 2, FITC-labeled)

Instruction Manual

Catalog Number	PK-AB704-60196F
Quantity	100 tests
Clone	2G11
IgG Subtype	mouse IgG1
Purity	> 95% IgG1 as determined by SDS-gel electrophoresis prior to addition of stabilizer and preservative.
Specificity	Specifically recognizes human p185HER-2.
Cross-reactivity	There are no data about cross-reactivity with other species.
Formulation	PBS (pH 7.4) with 0.1% (w/v) sodium azide
Storage	Stable at 2-8°C. After opening, aliquote contents and freeze at -20°C for long-term storage. Avoid repeated freeze/thaw cycles.
Shipping conditions	2-8°C
Applications	<p>Immunohistochemistry: PK-AB704-60196 can be used to stain cryostat sections or cell smears. Recommended for PK-AB704-60196 is the alkaline phosphatase-anti-alkaline-phosphatase (APAAP), or peroxidase anti-peroxidase (PAP) procedures or the three stage immunoperoxidase technique on acetone-fixed cryostat sections. PK-AB704-60196 may be used at a concentration of 1-5 µg/ml for cryostat sections.</p> <p>Flow cytometry: PK-AB704-60196 is also suitable as primary antibody in staining for FACS analysis, PK-AB704-60196F (FITC-labeled antibody) can be used for direct staining of cells, PK-AB704-60196B can be used in a two-step procedure using the biotin-(strept-)avidin system. Recommended dilution is 1:20 as determined on U937 cells (for exact concentration see vial label).</p> <p>Optimal antibody concentrations should be determined by the individual laboratory for each application.</p> <p>This antibody is also available as biotinylated conjugate.</p>

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.