

Data Sheet

p110α/p55γ

Human, Recombinant, Complex

Catalog #: 40645

Lot #: 161215 **Conc.:** 0.1 mg/ml

Formulation: 50 mM sodium phosphate, pH 7.0, 300 mM NaCl, 150 mM imidazole, 0.1 mM PMSF, 0.25 mM DTT, 25% glycerol.

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.

References:

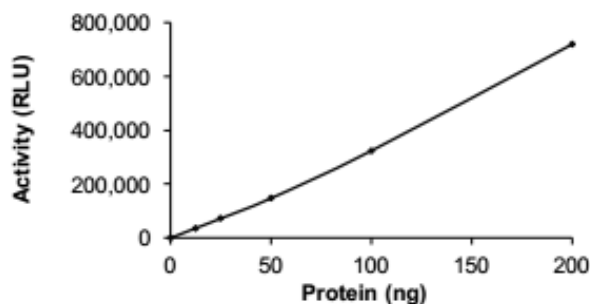
1. Terauchi, Y., *et al. Nature Genet.* **21:** 230-235 (1999)
2. Fukao, T., *et al., Nature Immunol.* **3:** 295-304 (2002)

Description: Complex of human p110α, also known as PIK3CA, GenBank Accession No. NM_006218, a.a. 2-1068(full length), with, N-terminal His-tag, MW=111 kDa, and human p55γ, GenBank Accession No. NM_003629.2, a.a. full length MW=61 kDa. Coexpressed in a Baculovirus infected Sf9 cell expression system.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

Specific Activity



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
p110α
and p55γ
Lane 2:
Protein Marker

MW: 111 kDa, p110α
61 kDa, p55γ
Purity: ≥70%

