

Data Sheet

p110α(H1047R)/p85α

Human, recombinant, FLAG-tag

Catalog #: 40641

Lot #: 131101-G **Conc.:** 0.2 mg/ml

Formulated in: 25 mM Tris, pH 8.0, 69 mM NaCl, 1.35 mM KCl, 0.03% Tween-20, 3 mM DTT, and 50% glycerol.

Stability: >6 months at -80°C. Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

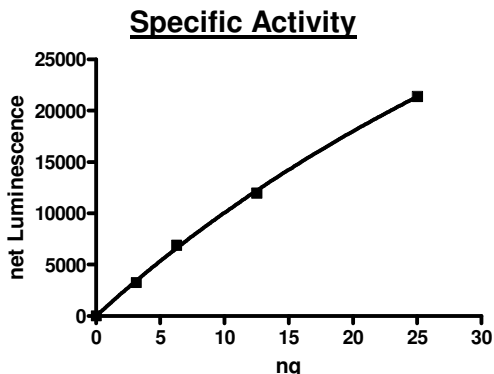
1. Engelman, J.A., *et al.* *Nature Med.* **14**: 1351-1356 (2008)
2. Hafner, C., *et al.*, *Proc. Natl. Acad. Sci. USA.* **104(33)**:13450-13454 (2007)
3. Serra, V., *et al.*, *Cancer Res.* **68(19)**:8022-8030 (2008)

Description: Complex of N-terminal FLAG-tagged recombinant full-length human p110α (GenBank Accession No. U79143), with H1047R mutation, and recombinant full length, human p85α (no tag, GenBank Accession No. XM_043865). Co-expressed in Sf9 cells in a Baculovirus expression system. p110α(H1047R) MW=125 kDa, p85α MW=84 kDa.

Specific Activity: 1770 pmol/min/μg. Assay Conditions: 25 μl kinase reaction is conducted in 40 mM Tris-HCl (pH 7.5), 10 mM MgCl₂, 100 μM ATP, 200 μM of PI3K substrate (BPS Cat.# 40560) and the PI3K enzyme for 18 minutes at 30°C. Conversion of ATP to ADP is detected using ADP-Glo[®] Luminescent Kinase Assay reagents (Promega Corp., Madison, WI).

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

Quality Assurance



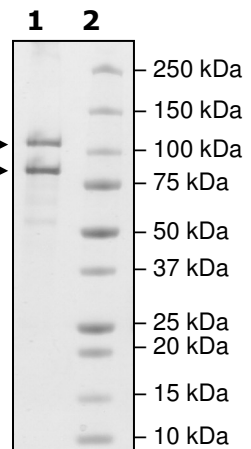
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
4.4 μg_p110α(H1047R)/p85α

Lane 2:
Protein Marker

MW:
p110α(H1047R):125 kDa
p85α: 84 kDa

Purity: ≥64%



OUR PRODUCTS ARE FOR RESEARCH ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
 To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com