

Product Information

| | |
|---------------------------|---|
| Construct: | KRAS (G12R) (His-2-185) |
| Mutation: | G12R |
| Concentration: | 2.41 mg/ml |
| Species: | Human |
| Formulated In: | 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 3 mM DTT. |
| Expression System: | <i>E. coli</i> |
| Format: | Aqueous buffer solution |
| Stability: | At least 6 months at -80°C. Avoid freeze/thaw cycles. |
| Storage: | -80°C |
| Genbank Accession: | NM_004985 |
| MW: | 23 kDa |
| Purity: | >90% |

Quality Control Data

4-20% SDS-Page Coomassie Staining

