

Data Sheet

JARID1B (PLU-1, KDM5B)

Mouse, recombinant, N-terminal FLAG-tag
Catalog #: 50122
Lot#: 100909 **Conc.:** 0.23 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 90 µg/ml, 10% glycerol, and 3 mM DTT.

Stability: >6 months at -80 °C

References:

1. Yamane K, *et al.* (2007). *Mol. Cell* **25 (6):** 801–12.
2. Barrett A, *et al.* (2007). *Int. J. Cancer* **121 (2):** 265–75.

Description:

Mouse JARID1B, also known as PLU-1 and KDM5B (GenBank Accession No. NM_152895) (amino-acids 2-751) with N-terminal FLAG-tag, MW=86 kDa, expressed in Sf9 cells via a baculovirus expression system.

Application:

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

Quality Assurance

10% SDS-PAGE Coomassie staining

Lane 1:
4 µg JARID1B
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 86 kDa
Purity: ≥50%

