

Data Sheet

Dicer, His-Tag

Human, Recombinant, N-terminal His-Tag

Catalog #: 79083

Lot #: 181026 **Conc.:** 0.17 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04%
Tween20, 20% glycerol, 40mM imidazole

Stability: At least 6 months at -80°C . Avoid
freeze/thaw cycles. Protein may be diluted to
 $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at
 -80°C .

References:

1. Hutvágner, G., *et al.* *Science*. 2001;
293(5531): 834-838.
2. Bernstein, E., *et al.*, *Nat. Genetics*.
2003; **35:** 215-217 (2003).

Description: Human Dicer, also known as
Dicer 1, and Ribonuclease III, GenBank
Accession No. NM_030621, a.a. 2-1992
(end), with a N-terminal His-Tag, expressed
in an Sf9 cell expression system. MW =221
kDa.

Quality Assurance

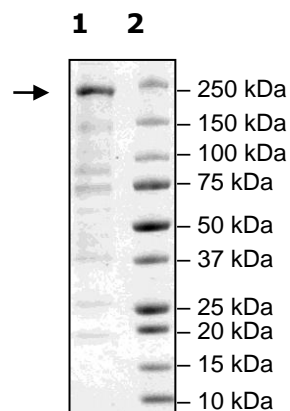
4-20% SDS-PAGE Coomassie staining

Lane 1:
Dicer
(2ug loaded)

Lane 2:
Protein Marker

Purity: $\geq 56\%$

MW:
221 kD



OUR PRODUCTS ARE FOR RESEARCH USE 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