

---

## Histone H3K79me2 Dimethyl Peptide

(Cat. No. R-1041)

---

|                                  |   |
|----------------------------------|---|
| <b>Description:</b>              | Synthesized dimethyl histone H3-K79 peptide derived from residues within 1-100 of human histone H3.   |
| <b>Purity:</b>                   | >90% by HPLC analysis   |
| <b>Concentration:</b>            | 0.5 mg/ml   |
| <b>Formulation:</b>              | Tris-buffered solution  |
| <b>Storage:</b>                  | -20°C, stable for 1 year from the date of shipment. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance. |
| <b>Application:</b>              | Enzyme activity<br>Histone affinity precipitation assay   |
| <b>Research use:</b>             | Research use only.  |
| <b>Handling Recommendations:</b> | For maximum recovery of the products, centrifuge the vial prior to opening the cap.   |

### Ordering Information

| Products         | Size   | Cat. No.   |
|------------------|--------|------------|
| Histone H3K79me2 | 100 µl | R-1041-100 |
| Dimethyl Peptide | 200 µl | R-1041-200 |

*This product is for research purposes only. Not intended for use in diagnostic procedures.*