

Cell lines & organisms analyzed with 5-EdU and 5-VdU (*de novo* DNA synthesis monitoring)

Cell line / Organism	Final 5-EdU concentration	Final 5-VdU concentration
HeLa cells	10 μM ^[1]	30 μM ^[11]
CEM cells ¹	20 μM ^[2]	
BT474 cells	0.1-20 μM ^[3]	
Jurkat cells	10 μM ^[4]	
NIH3T3 cells	10 μM ^[1]	
SK-BR-3 cells	0.1-10 μM ^[3]	
U2OS cells		30 μM ^[11]
MRC-5 cells		30 μM ^[11]
Vero cells		30 μM ^[11]
A549 cells		30 μM ^[11]
Primary human fibroblast (48BR (normal), 1BR (normal), XP15BR (XP-A), XP20BE (XP-G), XP13BR (XP-C), XP12BR (XP-D), CS10LO (CS-B))	10 μM ^[5]	
Mouse embryo	10-200 mg/kg ^[6]	
Chicken embryo	500 μM ^[7]	
Primary valvular interstitial cells	10 μM ^[8]	
Drosophila (neuroblasts, salivary glands, and wing discs)	15 μM ^[9]	
Plants (Alfalfa, Arabidopsis, grape, maize, rice and tobacco)	10 μM ^[10]	

Selected References:

- [1] Salic *et al.* (2008) A chemical method for fast and sensitive detection of DNA synthesis *in vivo*. *Proc. Natl. Acad. Sci. USA* **105**: 2415.
- [2] Hamelik *et al.* (2009) Click-iT™ Assay with Improved DNA Distribution Histograms. *Cytometry A* **75A**:862.
- [3] Diermeier-Daucher *et al.* (2009) Cell Type Specific Applicability of 5-Ethynyl-2'-deoxyuridine (EdU) for Dynamic Proliferation Assessment in Flow Cytometry. *Cytometry A* **75A**:535
- [4] Buck *et al.* (2008) Detection of S-Phase cell cycle progression using 5'-ethynyl-2'-deoxyuridine incorporation with click chemistry, an alternative to using 5'-bromo-2'-deoxyuridine. *Biotechniques* **44(7)**:927.
- [5] Limsirichaikul *et al.* (2009) A rapid non-radioactive technique for measurement of repair synthesis in primary human fibroblasts by incorporation of ethynyl deoxyuridine (EdU). *Nucleic Acids Res.* **37(4)**:e31.
- [6] Zeng *et al.* (2010) Evaluation of 5-ethynyl-2'-deoxyuridine staining as a sensitive and reliable method for studying cell proliferation in the adult nervous system. *Brain Research.* **1319**:21.

Cell lines & organisms analyzed with 5-EdU and 5-VdU (*de novo* DNA synthesis monitoring)

- [7] Warren *et al.* (2009) Chick embryo proliferation studies using EdU labeling. *Dev Dyn.* **238(4)**:944.
- [8] Monzack *et al.* (2012) A time course investigation of the statin paradox among valvular interstitial cell phenotypes. *American Journal of Physiology.* **303(7)**:H903.
- [9] Gouge *et al.* (2010) Detection of S Phase in multiple *Drosophila* tissues utilizing the EdU labeling technique. *Dros. Inf. Serv.* **93**: 213.
- [10] Kotogány *et al.* (2010) A rapid and robust assay for detection of S-phase cell cycle progression in plant cells and tissues by using ethynyl deoxyuridine. *Plant Methods* **6**: 5.
- [11] Rieder *et al.* (2014) Alkene-Tetrazine Ligation for Imaging Cellular DNA. *Angew. Chem. Int. Ed.* **53**: 9168.