

Cell lines & organisms analyzed with 5-EU (*de novo* RNA synthesis monitoring)

Cell line / Organism	Final 5-EU concentration
A549 cells ^[1]	0.2 mM
BEAS-2B cells ^[2]	1 mM
HEK 293T cells ^[3]	1 mM
LR7 cells ^[4]	0.25-4 mM
NIH3T3 cells ^[3]	1 mM
murine ES cell line R1 ^[5]	1 mM
Granyocytes ^[6] (isolated from PBMCs)	1 mM
Mouse ^[3] (intestine, kidney, liver, spleen)	2 mg/mouse
Plants ^[7] (<i>Vicia faba</i>)	1 mM

Selected References:

- [1] Xing *et al.* (2012) Foreign RNA Induces the Degradation of Mitochondrial Antiviral Signaling Protein (MAVS): The Role of Intracellular Antiviral Factors. *PLoS ONE* **7(9)**:e45136.
- [2] Abe *et al.* (2012) Analysis of interferon-beta mRNA stability control after poly(I:C) stimulation using RNA metabolic labeling by ethynyluridine. *Biochem Biophys Res Com* **428(1)**: 42.
- [3] Jao *et al.* (2008) Exploring RNA transcription and turnover *in vivo* by using click chemistry. *Proc. Natl. Acad. Sci. USA* **105 (41)**: 15779. [5] Xing *et al.* (2012) Foreign RNA Induces the Degradation of Mitochondrial Antiviral Signaling Protein (MAVS): The Role of Intracellular Antiviral Factors. *PLoS ONE* **7(9)**:e45136.
- [4] Hagemeyer *et al.* (2012) Visualizing Coronavirus RNA Synthesis in Time by Using Click Chemistry. *Journal of Virology* **86(10)**: 5808.
- [5] Vizlin-Hodzic *et al.* (2011) SAF-A Forms a Complex with BRG1 and Both Components Are Required for RNA Polymerase II Mediated Transcription. *PLoS ONE* **6(12)**:e28049.
- [6] Feng *et al.* (2013) Impairment of FOS mRNA Stabilization Following Translation Arrest in Granulocytes from Myelodysplastic Syndrome Patients. *PLoS ONE* **8(4)**:e61107.
- [7] Winnicki *et al.* (2012) Behavior of RNAs transcripts during nucleolus assembly and disassembly in *Vicia faba* root meristematic cells under normal conditions and after colchicine treatment. *Acta Physiol Plant* **34**: 1401.