



CRISPR-based depletion for smarter RNA-seq.

NEXTFLEX™ Cas9-gRNA Depletion Enzymes

The NEXTFLEX Cas9-gRNA Depletion Enzymes selectively eliminates abundant transcripts, freeing sequencing depth for biologically meaningful reads. Unlike other approaches, Cas9-gRNA depletion cleaves unwanted molecules at a dsDNA stage of your library prep workflow, reducing off-target effects and preserving transcript diversity even with low inputs or degraded material.

The NEXTFLEX Cas9-gRNA Depletion Enzymes are workflow-agnostic, compatible with bulk and single-cell RNA library prep for short or long read sequencing.

Key benefits:

- **Better.** Post-library depletion avoids manipulating input RNA, which can lead to biases or loss of material. Improved depletion efficiency.
- **Faster.** Ready to use. Workflow requires 1-minute hands-on time, and one-hour incubation. Straightforward implementation without complex steps.
- **Cheaper.** Improve outcome without significantly increasing the cost of library prep.

Product options

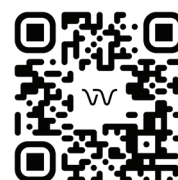
Target	Description	Species
rRNA	NEXTFLEX Cas9-gRNA Depletion Enzyme (Human/Mouse/Rat rRNA)	Human, Mouse, Rat
Globin	NEXTFLEX Cas9-gRNA Depletion Enzyme (Human Globin mRNA)	Human

Storage: -80°C | Shelf life: 12 months

How it works

The NEXTFLEX Cas9-gRNA Depletion Enzymes identify target DNA fragments within the NGS library and introduce precise double-strand breaks in the targeted RNA transcripts. Subsequent bead-based cleanup removes cleaved molecules, enriching for intact fragments prior to PCR amplification and sequencing. The result is cleaner libraries that yield more on-target reads and deeper coverage of transcripts of interest.

Transform your RNA-seq data with CRISPR-based depletion.
Learn more about the NEXTFLEX Cas9-gRNA Depletion Enzyme (Human/Mouse/Rat rRNA).



Learn more