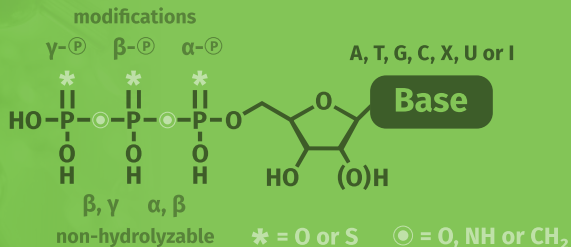


# NON-HYDROLYZABLE NUCLEOTIDES

Functional & Stable Analogs



# Non-hydrolyzable Nucleotides

## Functional & Stable Analogs

Discover our non-hydrolyzable nucleotide analogs (e.g., ATP $\gamma$ S, GppNHp)  
– ideal tools for studying signaling pathways, protein-nucleotide interactions and enzymatic activity without rapid hydrolysis.

### What we offer:

- More than 80 compounds ready to order as catalog products
- Custom synthesis option
- Direct contact with our scientists



Let's discuss  
your next project.