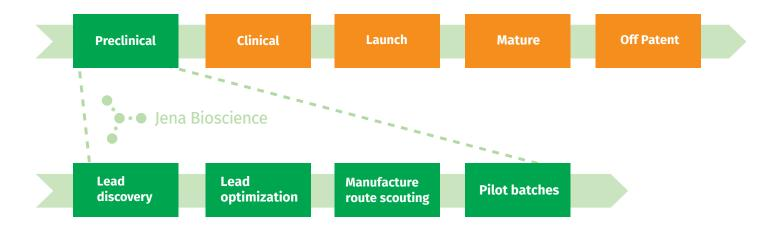


Jena Bioscience is a CRO active in chemical and enzymatic synthesis of nucleotides and their analogs. We have >1.500 compounds on stock, and our existing libraries of building blocks alongside with our specialized expertise allow short cut access to any desired novel nucleotide structure.



Outsourcing nucleotide chemistry projects to Jena Bioscience is an attractive option for clients from pharmaceutical and biotech industry as well as from academic research institutions.

## We offer

- > Very high success rates
- > Excellent track record
- > Speed and flexibility
- > Mile stone and / or FTE-based approaches from mg...kg scale
- > Attractive cost schemes affordable for academia

## Specialized nucleotide analogs

for pharmaceutical industry, biotech and academia





Jena Bioscience is a privately owned company founded in 1998 by a team of scientists from the Max-Planck-Institute for molecular Physiology in Dortmund. 30+ years of academic know how were condensed into the company in order to develop nucleotides and their analogs for medical and basic research.

Since the start up in 1998, the company has evolved into an established global reagent supplier with customers in 80+ countries. Our company premises are located in the city of Jena / Germany.

## Directly contact our people for more detailed information

## **Thomas Waldbach, PhD**Head of Chemistry Division Phone: +49-3641-628-5151

thomas.waldbach@jenabioscience.com



Sibylle Bauer, LL. M. oec.
Head of Marketing
Phone: +49-3641-628-5123
sibylle.bauer@jenabioscience.com

Thomas studied chemistry at the Friedrich-Schiller-University of Jena. During his graduate and postdoctoral research in Pretoria (South Africa) he specialized in small molecule organic synthesis which he continued at Roche before joining Jena Bioscience in 2001. Thomas was promoted Head of Chemistry in 2003 and has synthesized several thousands of different nucleotides during his career.

Sibylle studied law and economics at the Friedrich-Schiller-University of Jena and specialized in patent law of the biotech industry. After her bar exam she spent several internships in France before joining Jena Bioscience in 2003 as Head of Marketing.