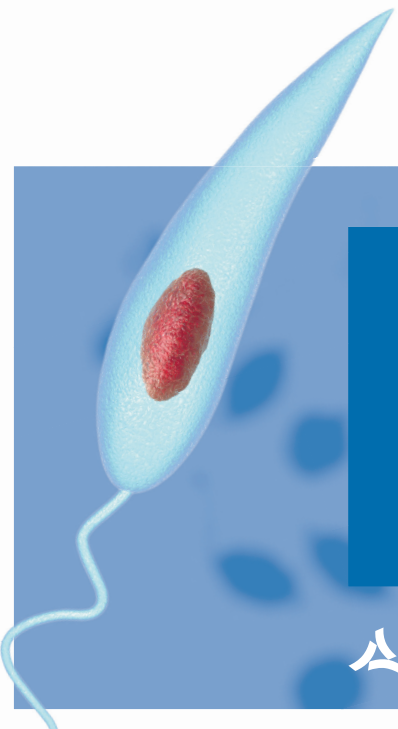




Jena Bioscience

Building Blocks of Life



The protozoan *Leishmania tarentolae* as a protein expression system

15.05.2018, 17 Uhr (HKI Kolloquium)

Dr. Andreas Licht (Jena Bioscience)

Hans-Knöll-Institut – Hörsaal Koch & Pasteur



IFTA AG
Certified QMS and EMS according to
DIN EN ISO 9001 and DIN EN ISO 14001
Reg.-No. ICV03597 034 and ICV03597 534

Lexsy Expression