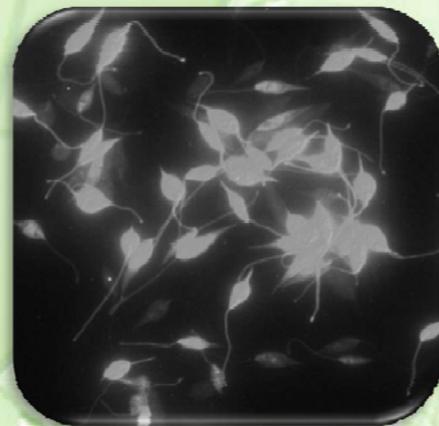




# LEXSY

## Eukaryotic Protein Expression with *Leishmania tarentolae*

- Biosafety Status -



Jena Bioscience GmbH  
Loebstedter Str. 80  
07749 Jena, Germany  
Tel.: +49-3641-628-5000  
Fax: +49-3641-628-5100

e-Mail: [info@jenabioscience.com](mailto:info@jenabioscience.com)



## LEXSY biosafety classifications

*Leishmania tarentolae* is classified as **biosafety level 1** in the USA.

Description on page 2.

Available via internet: Insert ATCC# „30143“ in

<http://www.lgcpromochem.com/atcc/>

*Leishmania tarentolae* is as well classified as S1

(Risikogruppe 1 = biosafety level 1) by the Central Biosafety Commission of Federal Republic of Germany (ZKBS).

Letter 10.12.1998 on page 3.

Mozilla

File Edit View Go Bookmarks Tools Window Help

Back Forward Reload Stop http://www.lgcpromochem.com/atcc/ Search

Home Bookmarks mozilla.org mozillaZine mozdev.org Jena Bioscience Google

**LGC Promochem**

About the ATCC-LGC Promochem Partnership

European Offices

How to Order

Special Forms

Technical Services

Products:

- Animal Virology
- Bacteria
- Bacteriophages
- Cell Lines & Hybridomas
- Molecular Genomics
- Fungi & Yeasts
- Plant Seeds
- Plant Virology
- Protistology
- Media, Reagents & Kits
- Special Collections
  - Yeast Genetic Research Resource Center
  - All Collections
  - Search Help
- View Your Shopping Cart

**Search**

Before submitting an order you will be asked to read and accept the terms and conditions of LGC's [Material Transfer Agreement](#).

**Protozoa And Algae**

**ATCC Number:** 30143

**Organism:** *Leishmania tarentolae* Parrot

**Designations:** Tar II

**Depositors:** SM Krassner

**Biosafety Level:** 1

**Isolation:** lizard, Tarentola mauritanica, Biakra, Algeria, 1934

**History:** ATCC<---SM Krassner <---W. Trager <--- L. Parrot

**Shipped:** frozen

**Propagation:** See Below

**Permits/Forms:** In addition to the [MTA](#) mentioned above, other ATCC and/or regulatory permits may be required for the transfer of this ATCC material. Anyone purchasing ATCC material is ultimately responsible for obtaining the permits. Please [click here](#) for information regarding the specific requirements for shipment to your location.

**Comments:** Mutant of W. Trager strain (ATCC 30267). Promastigotes. Effect of temperature on growth and nutritional requirements [4633] lipids [23748] Heterogeneity of minicircles [24027] isoenzyme electrophoresis for species identification [24036] Variation in kinetoplast minicircle and maxicircle DNA sequences [24219] Cyclopropane fatty acid [24043]

**Propagation:** ATCC medium: 1011 Diphasic blood agar medium  
Growth Conditions: axenic  
Temperature: 25.0 C  
See general instructions for thawing and storage of frozen material before proceeding. Add thawed contents to a single 10 x 125 mm glass screw-capped test tube of the appropriate medium. Incubate the culture vertically with the cap screwed on tightly. It is essential to establish cultures initially in small volumes. Once established, the culture can be scaled up to larger volumes. Vigorously agitate the culture and aseptically transfer 0.1 ml of culture to a fresh tube of medium weekly.

**References:** 4633: Krassner SM . Effect of temperature on growth and nutritional requirements of *Leishmania tarentolae* in a defined medium. J. Protozool. 12: 73-78, 1965. PubMed: 14292557  
23748: Beach DH , et al . Lipids of *Leishmania* promastigotes. J. Parasitol. 65: 201-216, 1979. PubMed: 448607

[Related Products](#)

