

Company Facts & Figures

Founded

1978

Founder

William A. Linton, Chairman and CEO

Governance

Independent and Privately Held

Number of Employees

1,400+ in 16 Countries

Number of Products

3,500+

Patent Leadership

Promega holds hundreds of patents in the areas of:

- Bioluminescence
- Cellular Analysis
- Cell biology
- Drug Discovery

Licensing

We offer licensing opportunities for a number of key technologies. The following technologies are available for licensing:

- NanoLuc® Luciferase
- NanoBIT® Technology
- HaloTag® Technology
- Green Fluorescent Protein
- Firefly Luciferase
- GloSensor™ Technology
- Apoptosis Marker Antibodies

Annual Revenue

Calendar Year 2016 Revenues Approaching 400M USD

Headquarters

Madison, WI, USA

Global Branches

Promega Corporation *Sydney, Australia*

Promega Biotecnologia do Brasil, Ltda. *São Paulo, Brazil*

Promega Biotech Co., Ltd. *Beijing, China*

Promega France *Lyon, France*

Promega GmbH *Mannheim, Germany (Austria/Poland)*

Promega Biotech India Private Ltd. *New Delhi, India*

Promega Italia, s.r.l. *Milan, Italy*

Promega KK *Tokyo, Japan*

Promega Korea, Ltd. *Seoul, South Korea*

Promega Benelux b.v. *Leiden, The Netherlands (Belgium/Luxembourg)*

Promega North America *Madison, Wisconsin, USA/Canada*

Promega Pte Ltd. *Singapore*

Promega Biotech Ibérica, SL *Madrid, Spain*

Promega Biotech AB *Stockholm, Sweden (Denmark/Estonia/Finland/Iceland/Norway)*

Promega AG *Dübendorf, Switzerland*

Promega UK *Southampton, United Kingdom*

Manufacturing

Promega Biosciences, LLC. *San Luis Obispo, CA, USA*

Promega BioSystems, Inc. *Seoul, South Korea*

Promega BioSystems Sunnyvale, Inc. *Sunnyvale, CA, USA*

Shanghai Promega Biological Products, Ltd. *Shanghai, China*

Over 50 Global Distributors