



42179949 Printed in USA. Rev. 11/17

Promega Campus Map

- ➊ **Agora.** Promega Headquarters. 5500 East Cheryl Parkway, Suite 110
- ➋ **Feynman Center.** Manufacturing. 2780 Woods Hollow Road
- ➌ **BioPharmaceutical Technology Center (BTC).** Manufacturing. 5445 East Cheryl Parkway
- ➍ **Research & Development Center (RDC).** 2800 Woods Hollow Road
- ➎ **Terso Solutions.** Subsidiary. 5540 Research Park Drive
- ➏ **Faraday Center.** Research & Development. 2800 South Fish Hatchery Road
- ➐ **Guest House & Conference Center.** 5474 Bjorksten Place
- ➑ **Earth House.** 5472 Bjorksten Place
- ➒ **da Vinci.** Maintenance & Engineering. 5511 Bjorksten Place

Additional Locations

- Aviation Operations.** Hangar. 3792 Corben Court, Madison, WI 53704
- Cell West.** Manufacturing. 8137 Forsythia Street, Middleton, WI 53562
- Kepler Center.** Operations. 5455 Nobel Drive, Fitchburg, WI 53711
- Rosalind Franklin Center (RFC).** Manufacturing. 2617 Progress Road, Madison, WI 53716
- 5500 Nobel Drive.** Instrumentation and Engineering. 5500 Nobel Drive, Fitchburg, WI 53711
- Mailing Address:** Promega Corporation. 2800 Woods Hollow Road, Madison, WI 53711-5399 United States. **Tel:** 608.274.4330. **Toll-Free:** 800.356.9526.