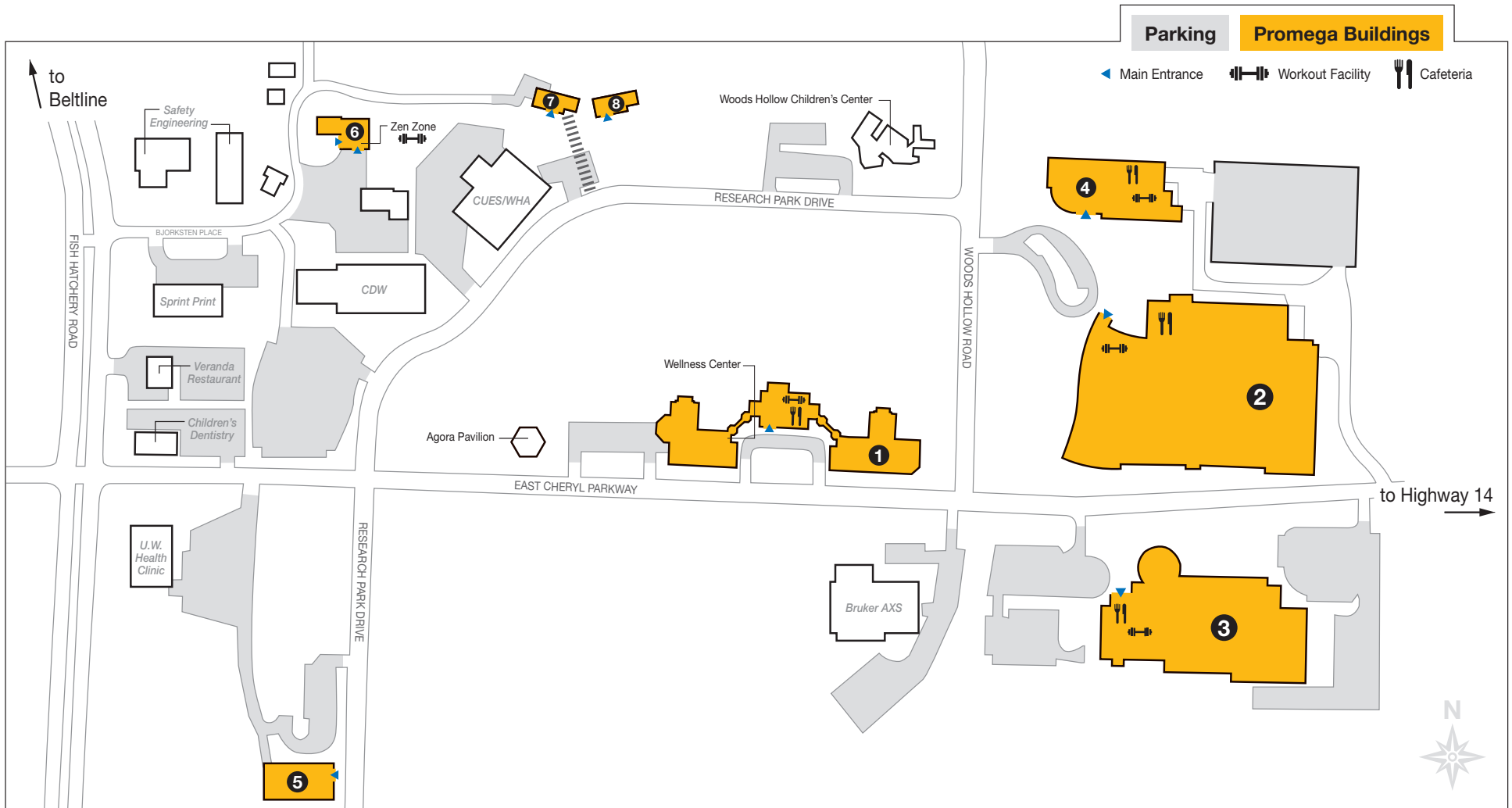




Promega Campus Map

Promega Corporation 2800 Woods Hollow Road Madison, WI 53711-5399 United States
tel: 608.274.4330 toll: 800.356.9526 www.promega.com



- 1** **Agora Promega Headquarters**
5500 East Cheryl Parkway, Suite 110
- 2** **Feynman Center Manufacturing**
2780 Woods Hollow Road
- 3** **BioPharmaceutical Technology Center (BTC) Manufacturing**
5445 East Cheryl Parkway
- 4** **Research & Development Center (RDC) R&D**
2800 Woods Hollow Road

- 5** **Terso Subsidiary**
5540 Research Parkway
- 6** **Faraday Center R&D**
2800 South Fish Hatchery Road
- 7** **Guest House & Conference Center**
5474 Bjorksten Place
- 8** **Earth House**
5472 Bjorksten Place

Not Shown:

- Rosalind Franklin Center (RFC) Production**
2617 Progress Road, Madison, WI 53716
- CPD Operations**
2920 Commerce Park Drive, Fitchburg, WI 53719
- Aviation Operations**
3792 Corben Court, Madison, WI 53704