



1180 E. Ellsworth Rd, Ann Arbor, MI 48108  
(800) 364-9897

### Directions to Cayman Chemical:

#### Detroit Metro Airport

Take I-94 W  
Take exit #177/State St  
Continue on S State St  
Turn left on Ellsworth Rd  
Arrive at 1180 E Ellsworth Rd on the Right

#### Ann Arbor Municipal Airport

Turn Left onto State Rd  
Turn Right onto Ellsworth  
Arrive at 1180 E Ellsworth Rd on the Right

#### Detroit Willow Run Airport

Turn Left on Airport Dr  
Turn Right on Tyler Rd  
Take ramp toward Wiard Rd  
Turn Right on Wiard Rd  
Take Left ramp onto Ecorse Rd [US-12]  
toward Ann Arbor  
Take ramp onto I-94 West  
Take exit #177/State St  
Continue on S State St  
Turn left on Ellsworth Rd  
Arrive at 1180 E Ellsworth Rd on the Right

