

1311 Orleans Drive  
Sunnyvale, CA  
94089-1136  
Ph: +1-800-635-5577  
Fx: +1-408-548-6461

## Hotel Information - Downingtown, PA

**Comfort Suites – Exton** - Corporate Rate *\$109.00*  
700 West Uwchlan Avenue, Exton, PA 19341, 610-594-4770  
[www.comfortsuites.com](http://www.comfortsuites.com)

- Complimentary breakfast served daily
- Wednesday evening managers reception - Beer/Wine/Soda/Appetizers
- Taxi Service to/from our office can be arranged by the front desk. When making your reservation please inform the reservation agent you will need taxi service.

**The Desmond Hotel** – Corporate Rate *\$157.00*  
One Liberty Blvd., Malvern, PA 19355, 610-296-9800  
[www.desmondgv.com](http://www.desmondgv.com)

- Hotel Shuttle Service to/from MDC Downingtown Office. When making your room reservation, please tell the reservation agent you will require hotel shuttle service and what time you will need it.

- 
- There is an Airport Shuttle Service to and from Philadelphia International (A2Z Airport Shuttle) for both hotels.  
<http://www.a2zairport.com/Schedule.asp>
  - If you prefer a town car, please call On Time Transport, 610-647-1354 (\$90.00 + gratuity – one way). They also provide taxi service as well.
  - If you choose to use a rental car, I recommend Enterprise (they have both an airport & Downingtown location), 1-800-261-7331.
  - Please note there is a Downingtown train station, however it is one mile from our office.