

The Alexandria at Torrey Pines

ADDRESS: 10996 Torreyana Road, San Diego, CA 92121



Parking & Driving Instructions

Parking Instructions

Parking is available to the general public off Callan Road.

Tenant Parking is available in the parking garage off Torreyana Road or available site parking.

To ensure the safety and proper use of our parking lot, please adhere to the following rules:

1. Remember to always lock your vehicle and remove any valuables, including cellular phones. Alexandria is not responsible for any damages to your vehicle, or theft.
 2. EV Charging is available to the public. Use is restricted to your time spent at The Alexandria. Please be considerate and move your vehicle out of EV Charging parking when not charging.
 3. Follow all marked parking space restrictions. ie. EV, ADA, etc.
 4. Notify Alexandria Concierge if you observe any hazards in the parking areas.
 5. Do not park illegally or in fire lanes.
-
1. Overnight parking is not permitted. Please notify the Alexandria Asset Services for extenuating circumstances.

2. Trailers and towed vehicles are not permitted in the parking areas.
3. Handicapped spaces are reserved for disabled persons only.
4. All vehicles must be parked in designated stalls and may occupy only one space.

**Cars parked found in violation to the parking rules are subject to citation and/or towing.*

Driving Directions

North going South on N Torrey Pines Road

- Going South on N Torrey Pines Road > follow right side ramp sign for Torrey Pines Golf Course The Lodge (ramp goes down) > turn left onto Callan Road and see the sign on the right for The Alexandria for outdoor parking > OR turn right onto Torreyana > turn right to the garage entrance
- *Public Transportation:* Please verify with Amtrak, Coaster and MTS schedule for availability

South going North on N Torrey Pines Road

- Going North on N Torrey Pines Road > turn R onto Science Park Road > turn L onto Torreyana Road > 3rd building on the left you will see a driveway to the front of the building parking or see the garage entrance > OR turn left on Callan Road and see sign on the left for The Alexandria
- *Public Transportation:* Please verify with MTS schedule for availability