

RESTAURANTS



SEA-GUINI @ OPAL SANDS \$\$\$
 430 South Gulfview Blvd. (727) 450-6236
 Waterview. Lunch 11am-3pm, Dinner 5pm-10pm.
 Beautiful views of the gulf from every seat, this Italian favorite features locally caught seafood, locally sourced produce and traditional pasta dishes all made from scratch in-house.
<http://www.seaguini.com/dinner/>



ISLAND WAY GRILL \$\$\$
 20 Island Way, Island Estates (1½miles) 727.461.6617
 Waterview. Mon-Thur 4pm-10pm, Fri & Sat till 11pm;
 Sat. Lunch 12Noon, Sunday Brunch 11-3pm, Dinner til 10pm
 A display case presents the day's catch, chops & deserts and chefs masterfully roll sushi, shuck oysters and grill over a wood fired hearth. Dine on Pan-Asian fare inside or on the Tiki deck over the water, either way you'll enjoy Clearwater's trendiest venue.
www.islandwaygrill.com/uploaded_files/islandwaymenu-2013-06.pdf



HEILMAN'S BEACHCOMBER \$\$\$
 447 Mandalay Avenue (½mile) 727.442.4144
 11:30am-10pm, Fri & Sat till 1030pm.
 Over 50 years of fine dining in one of the nicest dining rooms on Clearwater Beach, serving steaks, seafood & their signature Pan Fried Southern Chicken! Elegant, subdued setting popular with the classic crowd. Relax in plush comfort as the pianist plays your favorite tune.
www.heilmansbeachcomber.com/menus/dinner.html



COLUMBIA RESTAURANT \$\$\$
 1241 Gulf Blvd. Sand Key (2miles) 727.596.8400
 Waterview. 11:30am-10pm. The sister restaurant to the original 1905 in Ybor City. Feast on Spanish cuisine prepared using 100 year old recipes handed down through generations of the Hernandez & Gonzmart families. Enjoy casual dining and the panoramic view of the bay from indoors or out on the deck.
<http://www.columbiarestaurant.com/Menus-By-Location/Locations/Sand-Key.aspx>



CLEAR SKY CAFÉ \$\$
 490 Mandalay Avenue (¾mile) 727.442.3684
 7am-11pm, breakfast till 11:30am, Sat & Sun till 1pm.
 Breakfast, Lunch & Dinner. Happy hour 3-7pm daily, live acoustic guitar music on the patio nightly. Beach casual without being beach-bar rowdy. Great for families, wide-ranging menu has something for everyone.
www.clearskybeachsidecafe.com/Lunch-&-Dinner-Menu.html



CARETTA'S @ THE SANDPEARL RESORT \$\$\$\$
 500 Mandalay Avenue (¾mile) 727.674.4171
 Gulf-Front Waterview. Dinner served 5:30-10pm.
 Fine dining with a direct west-facing gulf view to enjoy the sun set. Fresh Florida cuisine, featuring a daily selection of local seafood, sushi bar, a wood burning oven, and a wine and cheese room. Extensive wine list with selections from around the world. Great for special occasions.
<http://www.sandpearl.com/dining/caretta-on-the-Gulf/>



FORLINI'S RISTORANTE \$\$\$
 435 Mandalay Avenue (½mile) 727.445.1155
 5-10pm, **CLOSED MONDAYS**
 Traditional family owned/operated white linen dining room, mornings will find mamma in the kitchen rolling fresh pasta & stuffing manicotti. Try the Zuppa di Pesce, Veal Saltimbocca or Filet Cacciatore and savor a port or cappuccino after at the dark wood bar. www.forlinis.com/menu_files/Menu.pdf



CESARE'S ITALIAN RESTAURANT \$\$\$
 794 South Gulfview Blvd. (¾mile) 727.449.8797
 Open 7days, 5-10pm.
 Classic pastas, pink sauces, meltingly soft gnocchi & a veal supper Frank Sinatra would sing for. Sample a home-made canoli or a crème brulee for desert. Dine inside or al-fresco, Gentlemen waiters provide professional, smart, attentive service.
www.cesareatthebeach.com



BOBBY'S BISTRO \$\$
 Behind Heilman's Beachcomber (½mile) 727.446.9463
 Open 7days, 5pm ...
 The less formal, less expensive "little brother" behind the Beachcomber. Great salads, pizza, pasta, fish, chops & daily specials. Dine in the open-air main dining room or the private leather booths in the Foxy Rock room. Say hello to the chef and see what's cooking as you walk in.
www.bobbysbistro.com/bistromenu.html



GUPPY'S (6miles) \$\$
 1701 Gulf Blvd., Indian Rocks Beach 727.593.2032
 11:30am-10pm, Fri. & Sat. till 10:30pm
 Where locals go for fresh fin & shell fish. Feast on fresh tuna, mussels, scallops or oysters - my favorite is the Peppercorn Tuna rare. Selections for landlubbers too. Family atmosphere in a converted cottage setting with tables inside or out on the front porch. Aquariums & consignment artwork add a colorful touch. Great for a light lunch.
www.3bestchefs.com/guppys/menus



E & E STAKEOUT (6miles) \$\$\$
 100 N. Indian Rocks Rd. Belleair Bluffs 727-585-6399
 Lunch Mon.-Fri. 11:30am-4pm, Dinner Sun.-Thur. 4pm-10pm, Fri. & Sat. till 10:30pm
 Sister restaurant to Guppy's, this steak house serves only the highest quality graded beef, along with options from the sea, in a zen-like atmosphere of rock, wood and water décor. Relax in plush leather semi circle booths or traditional tables and permit the staff to prepare a legendary feast. The spirit of Kokopelli lives!
www.3bestchefs.com/ee/menus



MARLIN DARLIN (6miles) \$\$
 2819 West Bay Drive, Belleair Bluffs, 727.584.1700
 Mon-Thur 4-10pm, Fri & Sat till 11pm, Sundays 11AM-10PM
 From the owners of the Island Way Grill & Salt Rock Grill, enjoy seafood and steaks in a dining space akin to a sport fishing yacht, with fighting chair bar stools and wood decked flooring. The stylish lounge and open-air kitchen resonate a vibrant atmosphere. Oysters, king crab & lobster tails, the Marlin Darlin is a prized catch.
www.marlindarlinkeywestgrill.com/menus.html



BONEFISH GRILL (6miles) \$\$
 2939 West Bay Drive, Belleair Bluffs, 727.518.1230
 4pm-10:30pm, Fri. & Sat. till 11:30pm, Sunday till 10pm.
 Zagat Rating 22. Part of the OSI Restaurant Group, featuring daily fresh fish specials prepared your choice of 4 different ways. Great appetizers, fresh fruit mojitos & martinis. Knowledgeable, attentive service. Not your average chain restaurant, here you're treated like family.
www.bonefishgrill.com/locator/details/belleair-bluffs-florida



SALT ROCK GRILL (10miles) \$\$\$
 19325 Gulf Blvd., Indian Shores, 727.593.7625
 Sun-Thur. 4pm-10pm, Fri. & Sat. till 11pm
 Zagat Rating 24, Wine Spectator Award, St.Pete Times Best Steak, Florida Beaches' Best Award winner. Hip atmosphere featuring monstrous steaks, crab & seafood. Pass a wood burning hearth, meat & fish display case and glass floor showcasing the wine cellar below on your way to the tiered dining area overlooking the water. Waterside bar & raw bar lounge offer a quiet respite.
www.saltrockgrill.com/menus.html



RUMBA ISLAND GRILL (5miles) \$\$
 1800 Gulf to Bay Blvd.(US 60), Clearwater 727.446.7027
 11:30am-9:30pm, Fri. & Sat. till 10pm, \$\$
 From the Island Way/Salt Rock Grill family of restaurants, come savor the jerk flavors of Jamaica and the islands of the Caribbean. Fresh fish with kinky aiolis, tropical salsas, big squeezes of lime or a splash of coconut milk. Island music, Island ambience - Yah Mon, Jamaican me hungry!
www.rumbaislandgrill.com/uploaded_files/rumba_menu-2014-06.pdf



BLACK PEARL (6miles) \$\$\$
 315 Main Street, Dunedin 727.734.3463
 Sun.-Thur. 5pm-11pm, Fri. & Sat. till midnight, \$\$\$
 Zagat Rating 26, St.Pete Times Best Splurge Night Out, Florida Beaches' Best Award winner. New American cuisine smartly prepared & expertly served in a cozy 12 table, intimate dining room reminiscent of SoHo Manhattan. Try the duck or filet, but save room for the Black Pearl ice cream. A real gem, perfect for that special occasion. Reservations required.
<http://www.theblackpearldunedin.com/menu/#everyday-menu>



CASA TINA (6miles) \$\$
 365 Main Street, Dunedin 727.734.9226
 11am-10pm, Fri. & Sat. till 11pm. \$\$
 Visit Tina's house and feast on fresh, healthy, authentic Mexican cuisine prepared by her husband Javier and served by her fabulous staff. Sauces, rice & beans and vegetarian dishes are prepared fresh daily. View art & artifacts they've collected on their travels through Mexico and enjoy a celebration of food & life. Visit Pan y Vino, their wine shop & living room next door, and relax after dinner over a game of backgammon or cribbage.
www.casatinas.com/menus.html



VILLA GALACE (7½miles) \$\$\$
 109 Gulf Blvd., Indian Rocks Beach 727.596.0200
 4:30pm-10:30pm. \$\$
 Prepare to become a member of 'The Family'. Receive a hearty handshake from Luigi as he greets you, seats you and treats you to some of his fresh-made pasta, or select from a menu featuring Mama's best recipes for seafood, veal, chicken and steak. Save room for cappuccino and desert. Professional service from dapperly dressed waiters.
www.villagallace.com/menu.html



GREEK TOWN GRILLE (3½miles) \$\$
 1222 Cleveland Street, Clearwater 727.447.7800
 Mon-Wed. 11:30am-10pm, Thur-Sat. 11:30am-2am, Sunday Noon-10pm. \$\$
 Lemony char-grilled octopus, crispy fried smelts with a zippy pepperoncini sauce, moussaka, pastitsio, chicken pita. My favorite are the lamb chops! Finish with a dessert called kantaifi, you will be dancing like Zorba afterwards. Great happy hour - Opa baby!
www.greektowngrille.com/menus

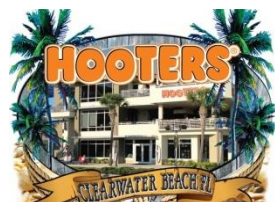
BEACH BARS



The Sand Bar @ Opal Sands

430 South Gulfview Blvd. 727.450.0380

Located a short walk down the dune off the south end of Clearwater Beach, Sandbar provides a relaxed atmosphere for lunch or dinner. We offer a delicious variety of Florida casual fare and some of the best drinks on the beach.



Hooters & The Lookout

381 Mandalay Ave.

(727) 443-7263

www.originalhooters.com/locations/tampa-bay/clearwater-beach

If you can't make it over to the original Hooters in Clearwater where the empire began 30 years ago, visit their newest location here on the beach overlooking the Clearwater Roundabout for your fix of chicken wings, curly fries and orange shorty-shorts. Hooters is on the 2nd floor and The Lookout is up on 3 with a fantastic view of Pier60 and the sunset. Over 30 TVs and live music on weekends up in the Lookout.

Toucan's

431 Mandalay Ave.

(727) 461-4500

Sports Bar with tons of TVs including 2 mega TVs on the north wall. Wide-ranging menu with all the popular bar offerings... try the Toucan Crunch Chicken Fingers or their Jumbo Wings. Three hi-top tables on the sidewalk allow for great people & traffic watching, live music or DJ on most nights. Open past 2am.



Jimmy's Crows Nest @ Pier House 60

101 Coronado Drive, 727.683.0002

www.pierhouse60.com/jimmys-crows-nest.html

Rooftop bar on 10th floor of the Pier House 60 Hotel, a great place to watch the sunset, with fantastic views of Gulf, Marina & the Roundabout. Projection TV shows your favorite ball game HUGE on the side of the building.



Reflections @ Opal Sands

430 South Gulfview Blvd 727.450.0380

Come join us pool side overlooking the beautiful Gulf of Mexico. Enjoy personalized service as our attendants bring handcrafted cocktails right to your lounge chair. **HOTEL GUESTS ONLY**



Brown Boxer

483 Mandalay Ave.

Pelican Plaza 727.441.6000

www.thebrownboxer.com

The other Sports Bar further up Mandalay Ave. Over 20 Flat Screen Hi-Def TVs with DirecTV Sports Pkgs, UFC & Boxing Pay-Per-Views. Live Music on weekends & full menu until 2am. Daily specials: Mondays' All-U-can eat Snow Crab; Tues Pint Night; Wed Ladies Night 2-4-1. Younger crowd hangs out until 3am.



Cork & Brew

524 Mandalay Ave.

727.483.5990

Craft Beer and upscale wine bar offers small plate tapas meals uniquely prepared using a searing hot volcanic slate! A scrumptious selection of desserts, great place to wind down on the cushy couches out on the sidewalk. Guitarist on weekends.

Frenchy's Rockaway Grill

7 Rockaway Street

727-446-4844

www.frenchyonline.com/locations/frenchys-rockaway

The flagship of Frenchy's four restaurant empire sits directly on the sand, one of the best spots to have a margarita & grouper sandwich while watching the sun set. Worth the wait for a table on the deck under the umbrellas. Live music nightly, satellite TV to entertain, bare feet & bikinis welcome.



Palm Pavilion

10 Bay Esplanade

727-446-2642

www.palmpavilion.com

An old favorite directly on the beach, enjoy a cool drink & lite bites or a grouper wrap while listening to the evening's guitarist. Check out the early photos of Clearwater Beach and join the applause when the sun sets over the water.





The Reef Steak and Seafood Co
 752 South Gulfview Blvd. 727.240.1182
<http://www.thereefsteakandseafood.com/>
 Excellent local bar, just down the street from our resort. full liquor bar, beer and wine, and delicious appetizers.



Frenchy's Salt Water Cafe
 419 Poinsettia Ave.
 727-461-6295
www.frenchyonline.com/locations/frenchys-saltwater-cafe
 2 blocks in from the beach, when Frenchy's Rockaway is 'slammed', head over here to beat the crowd for the same great menu but without the gorgeous gulf view.



Frenchy's South Beach Café
 351 South Gulfview Blvd
 727-441-9991
www.frenchyonline.com/locations/frenchys-south-beach-cafe
 On south Clearwater Beach overlooking the Beach Walk promenade. Stop in for a bite, cool drink & gentle misting after soaking up the hot sun out on the beach.



Cooters
 423 Poinsettia Ave.
 Clearwater Beach, FL 33767
 727-462-2668
www.cooters.com
 Looking for crab legs? This place is crawling with them, with Mon & Tues as All-U-Can-Eat Crab Legs! Local hang out with nightly dinner specials, entertainment, darts and sports. Plenty of TVs, during football season, wear a Pittsburgh Steelers T-shirt to receive a wink from the owner.



Crabby's
 333 South Gulfview Blvd.
 727.608.2065
www.crabbybills.com/Clwtr.html
 On the Beach Walk promenade just south of the Hyatt, enjoy sunsets from the second floor over the beach. Part of the Crabby Bill's family of restaurants, the menu features the old favorites as well as new creations.



Fusion Cigar Lounge
 483 Mandalay Ave.
 (727) 330-7649
www.fusioncigars.com
 The best cigar bar on the beach, Daisy Fuentes has been known to stop by. Huge walk-in humidor, comfy padded wicker couches outside in the courtyard; beer & wine and games of chess & dominoes to enjoy along with your smoke. Wine bar next door has guitarist on weekends.



Jimmy's Fish House & Iguana Bar
 Holiday Inn Gulfview
 521 South Gulfview Blvd.
 727.446.9720
www.jimmysfishhouse.net
 Hotel restaurant with a tiki deck and nice view overlooking Sand Key Inlet. Live music on weekends during high season.



Shepherd's Tiki Bar & Lounge
 601-619 South Gulfview Blvd.
 727.441.6875
www.shephards.com/entertainment/tiki-bar



The Wave Night Club @ Shepherd's
www.shephards.com/entertainment/night-club
 The area's largest hot and cold All-U-Can Eat buffet, open for breakfast, lunch and dinner. At dinner, pile your plate high with crab legs and slices from the carving station. Lots of choices: salads, sushi, fried chicken, salads, deserts. Eat in or outside over the inlet. Munchies available at the Tiki Bar. Entertainment at 3 different venues: Reggae on the tiki deck, pop inside the air conditioned lounge and techno-rock DJ in the Wave night club after 10pm Wed-Sun., cover charge. Great people watching!



Clearwater Wine Bar & Bistro
 483 Mandalay Ave.
 727.446.8805
www.clearwaterwinebar.com
 Wine bar with courtyard tables featuring over 40 wines by the glass, cheeses, sandwiches, salads, mini-pizzas, tacos & burritos. Happy hour 5-7pm nightly, guitarist on weekends; monthly wine tastings & 'Yappy Hour' for dog lovers. **CLOSED MONDAYS.**