

# Complete Genome Sequence and Methyome Analysis of *Beggiatoa leptomitiformis* strains D-401 and D-402.

Alexey Fomenkov<sup>a\*</sup>, Zhiyi Sun<sup>a</sup>, Tamas Vincze<sup>a</sup>, Maria Orlova<sup>b</sup>, Brian P. Anton<sup>a</sup>, Margarita Yu. Grabovich<sup>b</sup>, Richard J. Roberts<sup>a</sup>.

<sup>a</sup>New England Biolabs, Ipswich, MA, USA

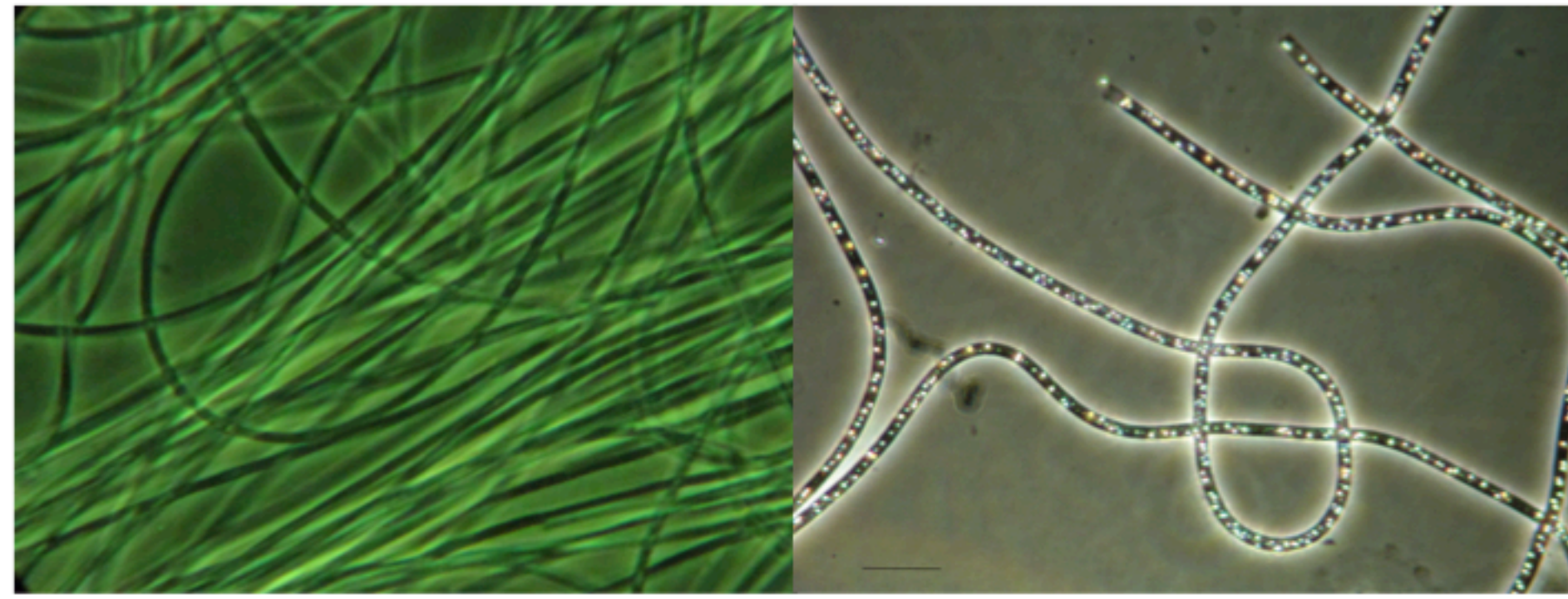
<sup>b</sup>Department of Biochemistry and Cell Physiology, Voronezh State University, Voronezh, Russia.

\*Correspondence: [fomenkov@neb.com](mailto:fomenkov@neb.com)

The taxonomy of *Beggiatoa* genus is still a work in progress. Despite many morphotypes of the *Beggiatoa* genus having been described in the literature, only one species, *B. alba* has been validated until now. In 2016, we described a second species, *B. leptomitiformis*. Two strains of *B. leptomitiformis* D-401 and D-402 have been isolated from different regions of Russia. They differ in their morphology and physiology and especially their ability to grow lithotrophically in the presence of thiosulfate. While *B. leptomitiformis* D-402 is able to accumulate elemental sulfur, strain D-401 cannot. We performed genomic sequencing of these two strains using the PacBio SMRT platform and assembled the reads into two complete circular genomes: *B. leptomitiformis* D-401 with 4,266,286 bp and D-402 with 4,265,296 bp. Both genome sequences have been deposited in GenBank with accession numbers CP018889 and CP012373 respectively (1). Surprisingly these two genomes showed almost 99% identity. The preliminary analysis of metabolic operons of thiosulfate oxidation (soxAXBZY) and autotrophic assimilation of CO<sub>2</sub> (the genes encoded for the enzymes of the Calvin-Bassham cycle) in both strains did not reveal any noticeable differences.

One advantage of the PacBio sequencing platform is its ability to detect the epigenetic state of the sequenced DNA, which allows for the identification of modified nucleotides and the corresponding motifs in which they occur. Thirteen DNA methyltransferase recognition motifs were found. They include one m4C and nine m6A modifications that were detected by direct SMRT sequencing and an additional three m5C motifs were detected in Tet2 treated DNA. The motifs were then matched with methyltransferase genes in the genome, and the results have been deposited in REBASE (2).

## Localization of Restriction and Modification genes on assembled chromosome map from *Beggiatoa leptomitiformis* strains.



The figure demonstrates micrograph of the bacterial filaments of *Beggiatoa leptomitiformis* strains D-401 without and D-402 with granules of elemental sulfur. The size of this bacterium can be as long as 100 micron and even longer. Bar on a right figure, 10um.

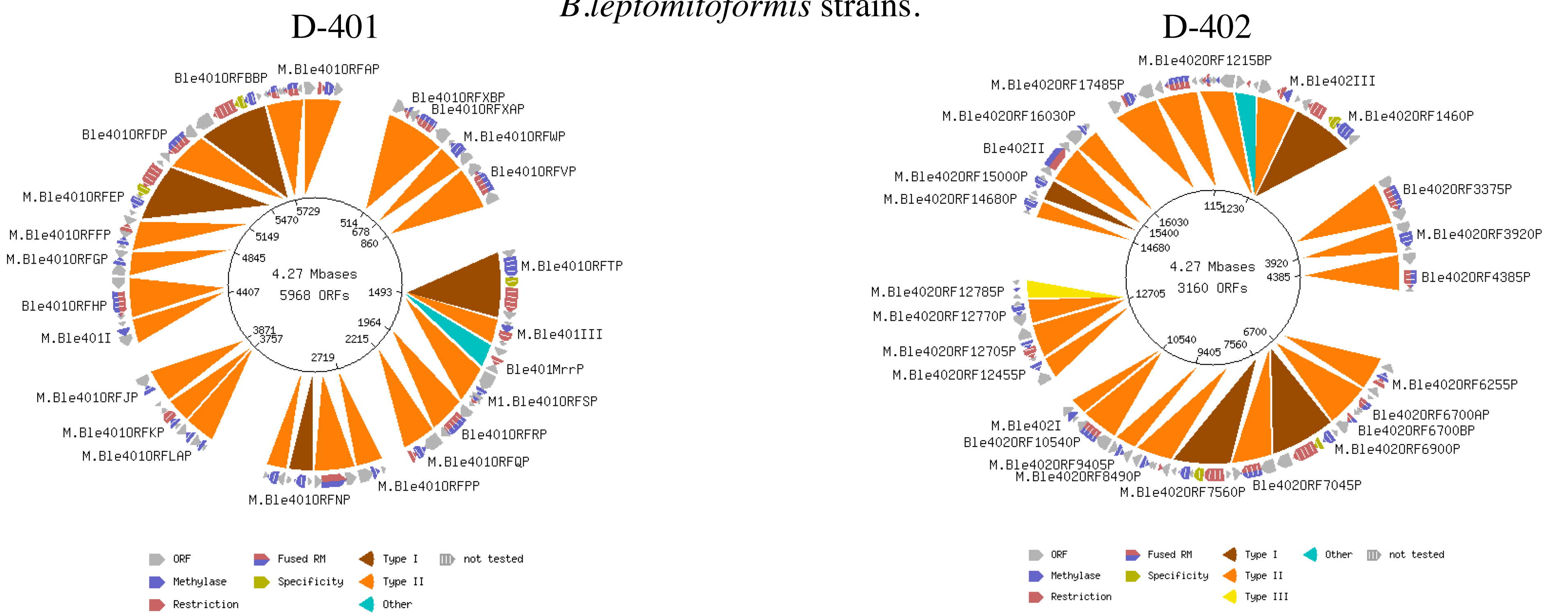


Table2 Methyltransferase genes and their cognate motifs identified in *B. leptomitiformis* strains D-401 and D-402

TYPE	LENGTH	COORDS	GENE AFFECTED	GENE COORDS	INTERNAL DELETION	HYPOTHETICAL PROTEIN
deletion	12064	12064	IAL038_RS00099	11777..12115	internal deletion	hypothetical protein
deletion	47013	47013	IAL038_RS00299	46477..47267	internal deletion	hypothetical protein
deletion	782	6181..6192	IAL038_RS00133	61269..62270	internal deletion	diguanylate cyclase response regulator
deletion	1441052	441052	IAL038_RS01885	439926..441278	internal deletion	magnesium transporter
deletion	719028	719028	IAL038_RS01329	719189..719664	internal deletion	hypothetical protein
deletion	782	89937..90018	IAL038_RS03848	898334..900254	internal deletion	long-chain fatty acid-CoA ligase
deletion	1926216	926216	IAL038_RS03945	925616..928076	internal deletion	l-deoxy-D-xylulose-5-phosphate synthase
deletion	1106556	106556	IAL038_RS04489	1060612..1064711	internal deletion	hypothetical protein
deletion	1209976	1209976	IAL038_RS05100	1209202..1210313	internal deletion	mexicin ABC transporter permease
deletion	853	123192..123944	IAL038_RS05105	1231830..1233006	internal deletion	hypothetical protein (arylsulfanyl nitrato)
deletion	782	147618..1477299	IAL038_RS06665	1476218..1476664	internal deletion	hypothetical protein (BleCSDN protein homolog, hypothetical protein)
deletion	782	147618..1477299	IAL038_RS06670	1476916..1477611	internal deletion	urea ABC transporter ATP-binding subunit (Ure)
deletion	853	180362..180474	IAL038_RS07529	1803331..1803648	internal deletion	hypothetical protein
deletion	853	180362..180474	IAL038_RS07529	1801901..1804767	internal deletion	hypothetical protein (Helicase P1a essential for oriC/DnaX-independent DNA replication)
deletion	1	1932313..1932373	IAL038_RS08079	1931320..1932756	internal deletion	arabinose channel protein
deletion	2044685	2044685	IAL038_RS08545	2044536..2044994	internal deletion	ribonuclease H1
deletion	782	2254406..2256294	IAL038_RS09799	2254368..2256066	internal deletion	hypothetical protein (Alkaline phosphatase (EC 3.1.3.1))
deletion	853	2636742..2637594	IAL038_RS11000	2636625..2636819	internal deletion	peptidase
deletion			INTERGENIC			hypothetical protein (Probable transmembrane protein)
deletion	853	2636742..2637594	IAL038_RS11005	2636991..2637894	internal deletion	hypothetical protein (Serine phosphatase Rbl1, regulator of sigma subunit)
deletion	853	2808212..2809064	IAL038_RS11790	2808941..2809562	internal deletion	hypothetical protein
deletion	1	3181477..3181477	IAL038_RS13330	3181223..3182407	internal deletion	hypothetical protein
deletion	1	3201097..3201097	IAL038_RS13410	3200947..3201966	internal deletion	hypothetical protein
deletion	1	3202925..3202925	IAL038_RS13411	3201970..3202925	internal deletion	hypothetical protein
deletion	1	3629468..3629468	IAL038_RS15340	3627872..3629689	internal deletion	hypothetical protein
deletion	1	3964261..3964261	IAL038_RS16620	3963872..3964276	internal deletion	transposase
deletion	1	19525..19525	INTERGENIC			
deletion	1	1168050..1168050	INTERGENIC			
deletion	1	1486221..1486221	INTERGENIC			
deletion	1	1729112..1729112	INTERGENIC			
deletion	1	2911338..2911338	INTERGENIC			
insertion	67850	67850	ORF_1878	67407..67826	internal insertion	Heat shock protein 60 family co-chaperone GroES
insertion	68002	68002	ORF_1883	67991..68360	internal insertion	Alkaline phosphatase (EC 3.1.3.1)
insertion	68006	68006	ORF_1880	68221..68749	internal insertion	hypothetical protein
insertion	68006	68006	ORF_1879	68477..68830	internal insertion	hypothetical protein
insertion	69038	69038	ORF_1877	68788..69315	internal insertion	unannoted protein product
insertion	69174	69174	ORF_1877	68788..69315	internal insertion	unannoted protein product
insertion	69038	69038	ORF_1878	68881..69267	internal insertion	unannoted protein product
insertion	102043	102043	ORF_1822	101586..102968	internal insertion	ATP-dependent DNA helicase RblB
insertion	1152026	1152026	ORF_1878	1152026..1152026	internal insertion	DNA repair protein (RecA)
insertion	120935	120935	ORF_1693	120219..120187	internal insertion	Ribonuclease P protein component (EC 3.1.26.5)
insertion	223992	223992	ORF_1670	223116..223474	internal insertion	Two-component hybrid sensor and regulator
insertion	227611	227611	ORF_1670	227582..228060	internal insertion	Methylene tetrahydrofolate synthetase dehydrogenase (EC 1.5.99.9)
insertion	1	262243..262243	ORF_1623	257822..262667	internal insertion	Adenylate cyclase (EC 4.6.1.1)
insertion	268884	268884	ORF_1616	265125..268886	internal insertion	hypothetical protein
insertion	272568	272568	ORF_1695	272568..272568	internal insertion	ATP-dependent DNA helicase UvrD/Pxa
insertion	853	343046..343898	ORF_1475	342817..344271	internal insertion	Membrane protein
insertion	853	343046..343898	ORF_1477	343781..344221	internal insertion	Membrane protein
insertion	1	344551..344551	ORF_1478	344549..344875	internal insertion	hypothetical protein
insertion	853	403645..404497	ORF_1404	401686..403647	internal insertion	sensory his histidine kinase/response regulator
insertion	853	403645..404497	ORF_1401	401814..404590	internal insertion	Ferrous siderophore transport system, periplasmic binding protein TonB
insertion	853	403645..404497	ORF_1402	401966..404590	internal insertion	Ferrous siderophore transport system, periplasmic binding protein TonB
insertion	853	403645..404497	ORF_1403	403975..404325	internal insertion	Ferrous siderophore transport system, periplasmic binding protein TonB
insertion	1	408019..408019	ORF_1388	407818..408887	internal insertion	Signal peptidase (EC 3.4.21.89)
insertion	1	520754..520754	ORF_1290	520754..520754	internal insertion	ATP-Oxidoreductase (1.8...)
insertion	1	720562..720562	ORF_939	720492..722214	internal insertion	hypothetical protein
insertion	1	726653..726653	ORF_937	723961..727632	internal insertion	Glyoxalase (EC 2.7.1.2)
insertion	1	729876..729876	ORF_938	729876..729876	internal insertion	Acyltransferase (F6-S-oxidoreductase)
insertion	1	744204..744204	ORF_915	743694..744556	internal insertion	Adenylate cyclase (EC 4.6.1.1)
insertion	1	813170..813170	ORF_828	812907..813349	internal insertion	hypothetical protein
insertion	1	898865..898865	ORF_879	897186..899043	internal insertion	Long-chain fatty acid-CoA ligase (EC 6.2.1.3)
insertion	1	1004515..1004515	ORF_872	1004299..1005012	internal insertion	Glyoxyl transferase, group 1
insertion	1	1004515..1004515	ORF_873	1004406..1004991	internal insertion	Glyoxyl transferase, group 1
insertion	1	1118880..1118880	ORF_4120	1118721..1120220	internal insertion	DNA polymerase III alpha subunit (EC 2.7.7.7)
insertion	1	1131646..1131646	ORF_4106	1131644..1132146	internal insertion	Protein DVU_0532 (HMC operon ORF 5)
insertion	1	1131646..1131646	ORF_4117	1131645..1131833	internal insertion	Protein DVU_0532 (HMC operon ORF 5)
insertion	1	1200500..1200500	ORF_324	1200478..1200526	internal insertion	ABC transporter, ATP-binding protein
insertion	1	1200550..1200550	ORF_325	1200469..1200521	internal insertion	ABC transporter, ATP-binding protein
insertion	853	1232316..1232406	ORF_287	1232320..1232631	internal insertion	diguanylate cyclase/phosphodiesterase (GGDEF & EAL domains) with PAS/PAC sensor(s)
insertion	853	1232316..1232406	ORF_288	1232320..1232631	internal insertion	diguanylate cyclase/phosphodiesterase (GGDEF & EAL domains) with PAS/PAC sensor(s)
insertion	1	1236071..1236071	ORF_148	1236071..1236071	internal insertion	hypothetical protein
insertion	1	1317119..1317119	ORF_148	1316927..1318904	internal insertion	hypothetical protein
insertion	1	1412842..1412842	ORF_54	1412447..1412851	internal insertion	Periplasmic septal ring factor with murcin hydrolase activity EnvC/YAP
insertion	1	1412842..1412842	ORF_53	1412858..1412858	internal insertion	Periplasmic septal ring factor with murcin hydrolase activity EnvC/YAP
insertion	1	1777115..1777115	ORF_5531	1775029..1777428	internal insertion	Multidomain transpeptidase/transglycosylase (EC 2.4.1.129) (EC 3.4...)
insertion	1	183874..183874	ORF_5432	1838342..1839412	internal insertion	Nitrogen regulation protein NtrB (EC 2.7.13.3)
insertion	1	1840221..1840221	ORF_5432	1840221..1840221	internal insertion	putative exported protein
insertion	1	1906363..1906363	ORF_5326	1906302..1907150	internal insertion	Hypothetical ATP-binding protein UPP0042, contains P-loop
insertion	1	209122..209122	ORF_5608	2090599..2096308	internal insertion	Periplasmic thioredoxin oxidoreductase DsbB, required for DsbA reoxidation
insertion	1	209122..209122	ORF_5609	2090599..2096308	internal insertion	Periplasmic thioredoxin oxidoreductase DsbB, required for DsbA reoxidation
insertion	1	2120882..2120882	ORF_5032	2120616..2120960	internal insertion	hypothetical protein
insertion	1	2286088..2286088	ORF_4803	2284152..2288318	internal insertion	conserved repeat domain
insertion	1	2287956..2287956	ORF_4803	2284152..2288318	internal insertion	conserved repeat domain
insertion	853	2359960..2359960	ORF_4711	2359921..2359977	internal insertion	hypothetical protein
insertion	853	2359960..2359960	ORF_4709	2360353..2361792	internal insertion	Sulfonamide modifying factor I precursor (C-alpha-formylglycine-generating enzyme 1)
insertion	1	2474762..2474762	ORF_4548	2474762..2474762	internal insertion	Quinolone biosynthesis QnrR-Radical SAM
insertion	1	2490818..2490818	ORF_4523	2489602..2491497	internal insertion	hypothetical protein
insertion	1	2607228..2607228	ORF_4537	2606789..2607781	internal insertion	hypothetical protein
insertion	1	2637171..2637171	ORF_4538	2636320..2640066	internal insertion	UDP-glucose 4-epimerase (EC 5.1.1.22)
insertion	853	2676217..2676217	ORF_4544	2676289..2677860	internal insertion	Glucosyl transferase G precursor
insertion	1	2724666..2724666	ORF_4183	2724366..2724748	internal insertion	hypothetical protein
insertion	1	2724666..2724666	ORF_4182	2724217..2724528	internal insertion	hypothetical protein
insertion	1	2773084..2773084	ORF_4120	2772610..2773461	internal insertion	Protein NCS-glyoxalase methyltransferase PmtC, methylates polypeptide chain release factors RFI and RF2
insertion	1	2773084..2773084	ORF_4121	2772908..2773237	internal insertion	Protein NCS-glyoxalase methyltransferase PmtC, methylates polypeptide chain release factors RFI and RF2
insertion	1	2927917..2927917	ORF_3640	2918187..2928472	internal insertion	Excisionase ABC subunit B
insertion	378	2979930..2980707	ORF_3640	2978242..2980254	internal insertion	hypothetical protein
insertion	378	2979930..2980707	ORF_3641	2980226..2980434	internal insertion	hypothetical protein
insertion	1	2989467..2989467	ORF_3640	2989467..2989467	internal insertion	High-affinity carbon uptake protein HtrHtrB
insertion	1	2994493..2994493	ORF_3646	2994431..2995222	internal insertion	Branched chain amino acid transport ATP-binding protein LAG (EC 3.1.4.1)
insertion	1	3131012..3131012	ORF_3648	3130760..3131932	internal insertion	COX6 COX13b
insertion	1	3161406..3161406	ORF_3603	3161006..3161809	internal insertion	Sensory histidine kinase BaeS
insertion	1	3161406..3161406	ORF_3604	3161164..3161553	internal insertion	Sensory histidine kinase BaeS
insertion	1	3178260..3178260	ORF_3572	3177741..3178868	internal insertion	diguanylate cyclase/phosphodiesterase (GGDEF & EAL domains) with PAS/PAC sensor(s)
insertion	1	3230972..3230972	ORF_3504	3230140..3231036	internal insertion	hypothetical protein
insertion	1	3301634..3301634	ORF_3386	3301448..3301660	internal insertion	SSU ribosomal protein S3p (S3p)
insertion	1	3301634..3301634	ORF_3387	3301448..3301660	internal insertion	SSU ribosomal protein S3p (S3p)
insertion	1	3301634..3301634	ORF_3385	3301613..3301959	internal insertion	SSU ribosomal protein S3p (S3p)
insertion	1	3467358..3467358	ORF_1611	3467170..3467367	internal insertion	dTPD-4-hydroxymethyl-5-pyrimidine (EC 5.1.3.13)
insertion	1	3470668..3470668	ORF_1611	3470668..3470668	internal insertion	methyltransferase, F8M family domain protein